

# 2019 Instrumentation and Consumable Catalog



# Resolvex® A200

Part Numbers:

**A200 96: 253-1160**

## Resolvex® A200 96

Resolvex A200 Standalone work station for automated sample preparation.

The compact benchtop offers the one stop solution for automating sample preparation utilizing creation of multiple work flows, and programmable dispensing of up to 11 solvents. Along with its innovative positive pressure system leading to clean samples, improving accuracy, throughput and enhancing the Life time of your analytical instrument.

The A200 comes with an easy to use touch screen interface allowing for easy set up of multiple work flows. In addition the light curtain safety feature will release gas pressure when manifold is activated to prevent any injuries.



# Resolvex<sup>®</sup> A100

Part Numbers:

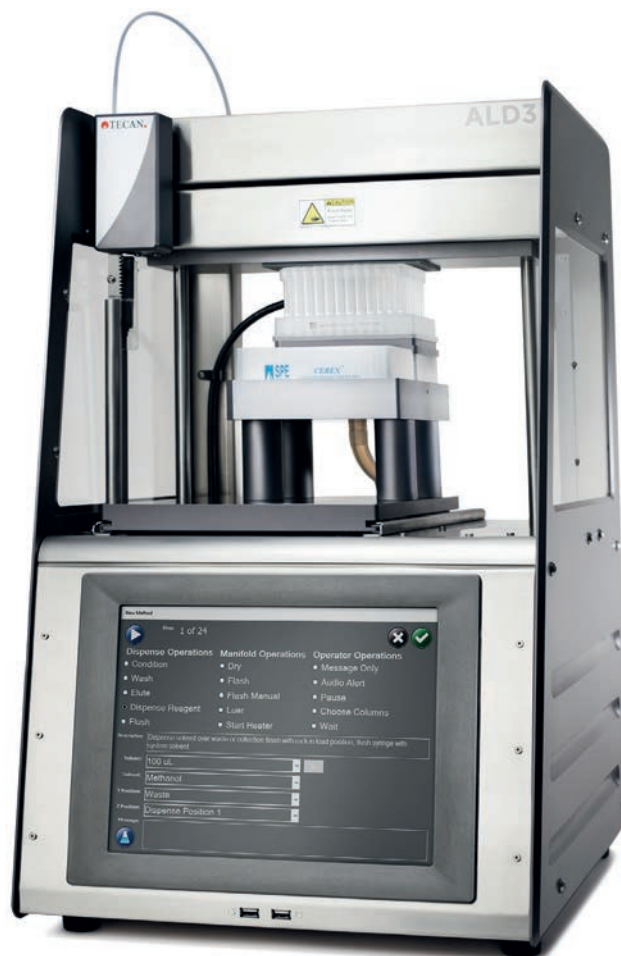
**A100 96: 253-0019**

**A100 48: 253-0014**

**Resolvex<sup>®</sup> A100 Standalone** work station for automated sample preparation.

The compact benchtop offers the one stop solution for automating sample preparation utilizing creation of multiple work flows, and programmable dispensing of up to 11 solvents. Along with its innovative positive pressure system leading to clean samples, improving accuracy, throughput and enhancing the Life time of your analytical instrument.

The A100 comes in 96 and 48 configuration allowing for for automated Work Flow solutions in multiple SPE formats.



# Resolvex® M10 96/M10 96 XT/M10 48

Part Numbers:

**M10 96 XT: 288-0006**

**M10 96: 288-0001**

**M10 48: 289-0004**

**Resolvex® M10 Standalone** work station for Positive Pressure solid phase extraction.

The manual Resolvex M10 48 and 96 are positive pressure manifold for 1, 3, and 6 ml cartridges, or 1ml 96 well plates. Restricted Flow gas ports allow single well pressure control for consistent processing of all samples.

The Resolvex M10 is simple to operate requires minimal training and can be deployed same day to help improving your work flow.

NOTE: The M10 96 comes in 2 configuration please consult your technical specialist to find the best configuration for your work.



# Cerex<sup>®</sup> C1 48 Concentrator 48

Part Numbers:  
**279-0050**

**Cerex<sup>®</sup> C48** is a 48 place evaporator that provides high throughput processing of sample extracts when used in conjunction with a 48 place A100 or Positive Pressure Processor.

Features:

- Clear Evaporation Chamber.
- Forced Air Heating - eliminates condensation problems from heated water baths.
- Smart Timer Control.
- Disposable Tips.
- 48 Individually Restricted Ports.
- Stepper Motor Driven - to adjust platform to any heights to different size tubes.





# Simplify your sample preparation.

## ADDITIONAL INSTRUMENTATIONS

Part Numbers	Descriptions
700-7001*	Evaporator Mini Vap STR. NEEDL. 110V
700-7004*	Evaporator Ultavap STR. NEEDLE 110V
282-0240	Column Heater 150 CERA
282-0252	Column Heater 250 CERA
282-0300	Column Cooler 250 CERA

\*Available in the US only

# Simplify your sample preparation.

## SPE COLUMNS PART NUMBER AND SORBENT DESCRIPTION



Material Sorbent Code	Sorbent Name	Example Applications
412	HP-SAX	THC/THCA
417	WWP2	Oasis HLB compatible sample preps
427	HP-SCX	Pain panal*, FNT, Aminoacids
606	PSAX-S	EtG/EtS, MMA
610	Trace-N	25-OH-VitD/steroids
611	Trace-N20	25-OH-VitD/steroids
613	Maestro	25-OH-VitD, Phospholipids Removal
615	PSAX	THC/THCA
633	OFX	Oral Fluid/Quantisal/Pain*
635	OFXQ	Oral Fluid/Quantisal/Pain*
637	OFQC	Pain/Oral-Eze
642	ETGS	EtG/EtS, MMA
667	FNT	Pain panal*, FNT, Aminoacids
671	SAX	Hair/THC/THCA
675	PWCX	Catecholamines
676	WWP	Plasma synthetic cannabinoids
682	THC	THC/THCA
691	CLN II	Pain panal*, FNT, Aminoacids
709	Trace-J	Pain panal*, FNT, Aminoacids
711	Trace-B	Pain panal*, FNT, Aminoacids
606/642	EtGS	MMA/EtG/EtS

**Pain Panel includes the following classes:**

Amphetamines, Barbiturates, Benzodiazepines, Carbamate, Cocaine, Cannabinoids, Gabapentinoids, Nicotine, Opioids, Phencyclidine, SSRIs, TCAs

**Amphetamines:**

Amphetamin, Methamphetamine, MDMA.

**Barbiturates:**

Allobarbital, Amobarbital, Aprobarbital, Alohenal, Barbital, Brallobarbital, Pentobarbital, Phenobarbitol, Secobarbitol, Thiopental.

**Phencyclidine:**

PCP, rolicyclidine (PCPy or 1-(1-phenylcyclohexyl)pyrrolidine); eticyclidine (PCE or N-ethyl-1-phenylcyclohexylamine); tenocyclidine (TCP or 1-(1-(2-thienyl)cyclohexyl)piperidine, 3-HO-PCP, 3-MeO-PCMo, and 3-MeO-PCP.

**SSRIs:**

Fluoxetine, Sertraline, Paroxetine, Escitalopram (Lexapro), Fluvoxamine, Citalopram, Volazodone (Viibrid), Vortioxetine, Seproxetine.

**TCAs:**

Butriptyline, Clomipramine, Imipramine, Trimipramine, Desipramine, Dibenzepin, Lofepamine, Maprotiline, Nortriptytine, Protriptyline, Amitriptyline, Amitriptylinoxide, Amoxapine, Demexiptiline, Dimetacrine, Dosulepin, Doxepin, Fluacizine, Imipraminoxide, Melitracen, Metapramine, Nitroxazepine, Noxiptiline, Pipofenzine, Quinupramine, Amineptine, Iprindole, Opipramol, Tianeptine.

**Opioids:**

Alfentanil, Buprenorphine, Carfentanil, Codeine, Dextropropoxyphene, Etorphine, Heroin, Hydromorphone, Methadone, Morphine, Naloxone, Norbuprenorphine, Oxycodone, Pethidine, Remifentanil, Sufentanil, Tramadol.

**Benzodiazepines:**

Alprazolam, Flunitrazepam, Chlordiazepoxide, Clonazepam, Diazepam, Lorazepam, Midazolam, Oxazepam, Prazepam, Lormetazepam, Nitrazepam, Temazepam, Zolpidam.

Tecan SP has not independently tested or verified performance of columns with all listed examples applications. Information presented is an aid to selection based on sorbent chemistry. Consistent with USP recommendations and good laboratory practice, the user must independently evaluate and validate: (a) the suitability of the method for their use, (b) their ability to process samples of their choosing following the method; and (c) their ability to proficiently perform the method in their facility with their personnel.



**HP-SAX**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polymeric Strong Anion Exchange	20 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
412-0021R-NBE	Column NBE HPSAX 1ml 2.5mg_96/pk
412-0051NBE	Plate HPSAX NBE 96 WELL 5MG _1/EA
412-0051NBEX	Plate HPSAX NBE 96 WELL 5MG _10PKS/CS
412-0101R	Column Cerex HPSAX 1ML 10MG _96/PK
412-0101RCH	Column Cerex HPSAX 1ML 10MG _1/EA
412-0201R	Column Cerex HPSAX 1ML 20MG _96/PK
412-0353C	Column Cerex HPSAX 3ML 35MG _100/PK
412-0353M	Column Cerex HPSAX 3ML 35MG _1000/PK
412-0401R	Column Cerex HPSAX 1ML 40MG _96/PK
412-0401RCH	Plate RCH CEREX HPSAX 1ML 96 40MG _1/EA
412-0401RCHX	Plate RCH HPSAX 1ML 96 40MG _10PKS/CS

**WWP2**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polymeric Reverse Phase	30 $\mu$	Water Wettable Polymer

<b>Part Numbers</b>	<b>Descriptions</b>
417-0051R-NBE	Column NBE WWP2 1ML 5.0MG _96/PK
417-0101R	Column Cerex WWP2 1ML 10MG _96/PK
417-0101RCH	Plate RCH Cerex WWP2 1ML 10MG _1/EA
417-0101RM	Column Cerex WWP2 1ML 10MG _960/PK
30147136	Column Cerex WWP2 3ML 35MG _100/PK
30147142	Column Cerex WWP2 3ML 35mg 1000/PK

**HP-SCX**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polymeric Strong Cation Exchange	30 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
427-0021NBEX	Plate NBE HPSCX 96 WELL 2.5MG _10PKS/CS
427-0021RM-NBE	Column NBE HPSCX 1ML 2.5MG _960/PK
427-0021R-NBE	Column NBE HPSCX 1ML 2.5MG _96/PK
427-0051NBE	Plate NBE HPSCX 96 WELL 5.0MG _1/EA
427-0051NBEX	Plate NBE HPSCX 96 WELL 5.0MG _10PKS/CS
427-0051RM-NBE	Column NBE HPSCX 1ML 5.0MG _960/PK
427-0051R-NBE	Column NBE HPSCX 1ML 5MG _96/PK
427-0101R	Column Cerex HPSCX 1ML 10MG _96/PK
427-0353C	Column Cerex HPSCX 3ML 35MG _100/PK
427-0353M	Column Cerex HPSCX 3ML 35MG _1000/PK
427-0506C	Column Cerex HPSCX 6ML 50MG _100/PK
427-0506M	Column Cerex HPSCX 6ML 50MG _1000/PK

**HP-PSAXS**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Strong Anion Exchange	10 $\mu$	Micro Particulate Silica

<b>Part Numbers</b>	<b>Descriptions</b>
606-0501R	Column C18 NON-EC 3ML 100MG _50/PK
606-0501RCH	Column Cerex PSAXS 1ML 50MG _96/PK
606-0503C	Plate RCH PSAXS 1ML 96 50MG _1/EA
606-0503M	Column Cerex PSAXS 3ML 50MG _100/PK

**TRACE N**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
C18 Endcapped	10 $\mu$	Micro Particulate Silica

<b>Part Numbers</b>	<b>Descriptions</b>
610-0101R	Column Cerex TRACE-N 1ML 10MG _96/PK
610-0101RCH	Plate RCH TRACE-N 1ML 96 10MG _1/EA
610-0101RCHX	Plate RCH TRACE-N 1ML 96 10MG _10PKS/CS
610-0101RM	Column Cerex TRACE-N 1ML 10MG _960/PK
610-0201R	Column Cerex TRACE-N 1ML 20MG _96/PK
610-0201RM	Column Cerex TRACE-N 1ML 20MG _960/PK
610-315C	Column Cerex TRACE-N 3ML 15MG _100/PK
610-315M	Column Cerex TRACE-N 3ML 15MG _1000/PK

**TRACE N-20**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
C18 Endcapped	15 $\mu$	Micro Particulate Silica

<b>Part Numbers</b>	<b>Descriptions</b>
611-0051NBE	Plate TRACE-N20 96 WELL NBE 5MG _1/EA
611-0051R-NBE	Column NBE TRACE-N20 5.0MG _96/PK
611-0101R	Column Cerex TRACE-N20 1ML 10MG _96/PK
611-0101RCH	Plate RCH TRACE-N20 1ML 96 10MG _1/EA
611-0101RCHX	Plate RCH TRACE-N20 1ML 96 10MG _10PKS/CS
611-315M	Column Cerex TRACE-N20 3ML 15MG _1000/PK

**MAESTRO**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
C18 Un-encapped	10 $\mu$	Micro Particulate Silica

<b>Part Numbers</b>	<b>Descriptions</b>
613-0051NBE	Plate NBE Maestro-A 96 WELL 5MG_1/EA
613-0051R-NBE	Column NBE Cerex Maestro A 1ML 5MG_96/PK
613-0151R	Column Cerex Maestro 1ML 15MG _96/PK
613-0151RCH	Plate RCH Cerex Maestro 1ML 96 15MG_1/EA
613-0151RCHX	Plate RCH Maestro 1ML 96 15MG _10PKS/CS
613-0151RM	Column Cerex Maestro 1ML 15MG _960/PK

**PSAX**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Strong Anion Exchange	15 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
615-0051NBE	Plate NBE PSAX 96 WELL 5.0MG _1/EA
615-0051NBEX	Plate NBE PSAX 96 WELL 5.0MG _10PKS/CS
615-0051RM-NBE	Column NBE PSAX 1ML 5.0MG _960/PK
615-0101R	Column Cerex PSAX 1ML 10MG _96/PK
615-0101RCH	Plate RCH Cerex PSAX 1ML 96 10MG _1/EA
615-0101RCHX	Plate RCH Cerex PSAX 1ML 10MG _10PKS/CS
615-0101RM	Column Cerex PSAX 1ML 10MG _960/PK
615-0201R	Column Cerex PSAX 1ML 20MG _96/PK
615-0201RCH	Plate RCH Cerex PSAX 1ML 20MG _1/EA
615-0353C	Column Cerex PSAX 3ML 35MG _100/PK
615-0353M	Column Cerex PSAX 3ML 35MG _1000/PK

**OFX**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Proprietary Ion Exchange	30 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
633-0031NBE	Plate NBE OFX 96 WELL 3MG _1/EA
633-0031NBEX	Plate NBE OFX 96 WELL 3MG _10PKS/CS
633-0031RM-NBE	Column NBE OFX 1ML 3MG _960/PK
633-0031R-NBE	Column NBE OFX 1ML 3MG _96/PK
633-0101NBE	Plate OFX 96-WELL NBE 10MG
633-0101R	Column Cerex OFX 1ML 10MG _96/PK
633-0353C	Column Cerex OFX 3ML 35MG _100/PK

**OFX - Q**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Proprietary Ion Exchange	30 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
635-0031NBE	Plate NBE OFXQ 96 WELL 3MG _1/EA
635-0031NBEX	Plate NBE OFXQ 96 WELL 3MG _10PKS/CS
635-0031RM-NBE	Column NBE OFXQ 1ML 3MG _960/PK
635-0031R-NBE	Column NBE OFXQ 1ML 3MG _96/PK
635-0051NBE	Plate NBE OFXQ 96 WELL 5MG _1/EA
635-0051NBEX	Plate NBE OFXQ 96 WELL 5MG _10PKS/CS
635-0051RM-NBE	Column NBE OFXQ 1ML 5MG _960/PK
635-0051R-NBE	Column NBE OFXQ 1ML 5MG _96/PK
635-0051R-NBE-HC	Column NBE OFXQ HC F. 1ML 5MG _96/PK
635-0101R	Column Cerex OFXQ 1ML 10MG _96/PK
635-0051R-NBE-HC	Column NBE OFXQ HC F. 1ML 5MG _96/PK
635-0101RCH	Plate RCH CEREX OFXQ 1ML 96 10MG _1/EA
635-0101RCHX	Plate RCH OFXQ 1ML 96 10MG _10PKS/CS
635-0101RM	Column Cerex OFXQ 1ML 10MG _960/PK

**OFQC**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Proprietary Ion Exchange	30 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
637-0051NBE	Plate NBE OFQC 96 WELL 5MG _1/EA
637-0051NBE-HC	Plate NBE OFQC 96WELL HC F. 5MG _1/EA
637-0051NBEX	Plate NBE OFQC 96 WELL 5.0MG _10PKS/CS
637-0051NBEX-HC	Plate NBE OFQC 96WELL HC F 5MG _10PKS/CS
637-0051RM-NBE	Column Cerex OFQC 1ML NBE 5MG _960/PK
637-0051R-NBE	Column NBE OFQC 1ML 5MG _96/PK
637-0101R	Column Cerex OFQC 1ML 10MG _96/PK
637-0101RCH	Plate RCH CEREX OFQC 1ML 96 10MG _1/EA

**ETGS**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Strong Anion	10 $\mu$	Micro Particulate Silica

<b>Part Numbers</b>	<b>Descriptions</b>
642-0501R	Column Cerex ETGS 1ML 50MG _96/PK
642-0501RCH	Plate RCH Cerex ETGS 1ML 96 50MG _1/EA
642-0501RCHX	Plate RCH ETGS 1ML 96 50MG _10PKS/CS
642-0501RM	Plate NBE OFQC 96WELL HC F 5MG _10PKS/CS
642-0503C	Column Cerex ETGS 3ML 50MG _100/PK
642-0503M	Column Cerex ETGS 3ML 50MG _1000/PK

**PWCX**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Strong Cation Exchange	15 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
675-0051R-NBE	Column NBE Cerex PWCX 1ML 5MG _96/PK
675-0201R	Column Cerex PWCX 1ML 20MG _96/PK
675-0201RCH	Plate RCH PWCX 1ML 96 20MG _1/EA
675-0351R	Column Cerex PWCX 1ML 35MG _96/PK
675-0351RCH	Plate RCH Cerex PWCX 1ML 96 35MG _1/EA
675-0353M	Column Cerex PWCX 3ML 35MG _1000/PK

**WWP**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polymeric Reverse Phase	15 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
676-0101R	Column Cerex WWP 1ML 10MG _96/PK
676-0101RCH	Plate RCH Cerex WWP 1ML 96 10MG _1/EA
676-0151RCH	Plate RCH Cerex WWP 1ML 96 15MG _1/EA
676-0201R	Column Cerex WWP 1ML 20MG _96/PK
676-0353C	Column Cerex WWP 3ML 35MG _100/PK
676-0353M	Column Cerex WWP 3ML 35MG _1000/PK

**THC**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polymeric Reverse Phase	15 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
682-0101R	Column Cerex THC 1ML 10MG _96/PK
682-0353	Column Cerex THC 3ML 35MG _500/PK
682-0353C	Column Cerex THC 3ML 35MG _100/PK
682-0353M	Column Cerex THC 3ML 35MG _1000/PK
682-0503	Column Cerex THC 3ML 50MG _500/PK
682-0503C	Column Cerex THC 3ML 50MG _100/PK
682-0506	Column Cerex THC 6ML 65MG _500/PK
682-0506C	Column Cerex THC 6ML 65MG _100/PK
682-0506M	Column Cerex THC 6ML 65MG _1000/PK



**PSCX**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polymeric Strong Cation Exchange	15 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
687-0021NBE	Plate NBE PSCX 96 WELL 2.5MG _1/EA
687-0021NBEX	Plate NBE PSCX 96 WELL 2.5MG _10PKS/CS
687-0021RM-NBE	Column NBE PSCX 1ML 2.5MG _960/PK
687-0021R-NBE	Column PSCX 1ML NBE 2.5MG _96/PK
687-0051NBE	Plate NBE Cerex PSCX 96 WELL 5MG _1/EA
687-0051NBEX	Plate PSCX NBE 96 WELL 5MG _10PKS/CS
687-0051RM-NBE	Column NBE PSCX 1ML 5MG _960/PK
687-0051R-NBE	Column PSCX NBE 1/EA 1ML 5MG _96/PK
687-0101R	Column Cerex PSCX 1ML 10MG _96/PK
687-0101RCH	Plate RCH Cerex PSCX 1ML 96 10MG _1/EA
687-0101RCH2	Plate RCH2 PSCX 1ML 96 10MG _1/EA
687-0101RCH2X	Plate RCH2 PSCX 1ML 96 10MG _10PKS/CS
687-0101RCH-48A	Plate RCH Cerex PSCX 48 1ML 10MG _2/PK
687-0101RCH-48C	Plate RCH CKRD PSCX 1ML 48 10MG _1/EA
687-0101RCHX	Plate RCH PSCX 1ML 96 10MG _10PKS/CS
687-0101RCHX-48A	Plate RCH Cerex PSCX 48 1ML 10MG _20/PK
687-0101RCHX-48C	Plate RCH CKRD PSCX 1ML 48 10MG _10PKS/CS
687-0101RM	Column Cerex PSCX 1ML 10MG _960/PK
687-0201R	Column Cerex PSCX 1ML 20MG _96/PK
687-0201RCH	Plate RCH Cerex PSCX 1ML 96 20MG _1/EA

**CLIN II**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polymeric Strong Cation Exchange	15 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
691-0353	Column Cerex CLIN II 3ML 35MG _500/PK
691-0353C	Column Cerex CLIN II 3ML 35MG _100/PK
691-0353D-HC	Column Cerex CLIN II HC F. 3ML _500/PK
691-0353M	Column Cerex CLIN II 3ML 35MG _1000/PK
691-0506	Column Cerex CLIN II 6ML 50MG _500/PK
691-0506C	Column Cerex CLIN II 6ML 50MG _100/PK
691-0506M	Column Cerex CLIN II 6ML 50MG _1000/PK

**TRACE - J**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polymeric Strong Cation Exchange	15 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
709-0101RCH	Plate RCH Cerex TRACE-J 96 1ML 10MG _1/EA
709-0101RCHX	Plate RCH Trace-J 1ML 96 10MG _10PKS/CS
709-0101RM	Column Cerex Trace-J 1ML 10MG _960/PK
709-0506C	Column Cerex Trace-J 6ML 50MG _100/PK
709-0506M	Column Cerex Trace-J 6ML 50MG _1000/PK
709-335C	Column Cerex Trace-J 3ML 35MG _100/PK
709-335M	Column Cerex Trace-J 3ML 35MG _1000/PK

**TRACE - B**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polymeric Strong Cation Exchange	15 $\mu$	PSDVB

<b>Part Numbers</b>	<b>Descriptions</b>
711-0101R	Column Cerex Trace-B 1ML 10MG _96/PK
711-0101RCH	Plate Cerex Trace-B 1ML 10MG _1/EA
711-0101RCHX	Plate Cerex Trace-B 1ML 10MG_10/CS
711-0101RM	Column Cerex Trace-B 1ML 10MG _960/PK
711-0201R	Column Cerex Trace-B 1ML 20MG _96/PK
711-0201RM	Column Cerex Trace-B 1ML 20MG _960/PK
711-0506C	Column Cerex Trace-B 6ML 50MG _100/PK
711-0506M	Column Cerex Trace-B 6ML 50MG _1000/PK
711-335C	Column Cerex Trace-B 3ML 35MG _100/PK
711-335M	Column Cerex Trace-B 3ML 35MG _1000/PK

**CFP**

<b>Sorbent Type</b>	<b>Particle Size</b>	<b>Base Material</b>
Polypropylene or PTFE	N/A	N/A

<b>Part Numbers</b>	<b>Descriptions</b>
961-0001CFP	Plate NBE CFP 96 WELL _1/EA (Polypropylene)
961-0001R-CFP	Column CFP 1 NBE _96/PK (Polypropylene)
962-0001CFP	Plate ND CFP 96 WELL NBE _1/EA (PTFE)
962-0001R-CFP	Column ND CFP NBE _96/PK (PTFE)

# Simplify your sample preparation.

## ACCESSORIES



### 96 WELL SPACER REPLACEMENT PART NUMBERS

Part Number	Description	Description 2	Guarantee Compatible Collection Plates*	Packaging
253-0038-SP	Waste Barriers low profile insert	Used with Waste Spacer	N/A	5/pk
253-0037-SP	Waste Barriers low profile insert	Aluminum spacer used with Waster barrier	N/A	Ea.
253-0023-SP	RCB-N1	Aluminum Support for the Cerex 96 well RCH products.	Nunc	Ea.
253-0105	NBE-N1 Spacer	Aluminum For NBE 96 Well	Nunc	Ea.
30137227	RCB-G1	Aluminum support for the Cerex 96 well RCH products	Nunc	Ea.

\*Note: Spacers and rack can also fit other brand of collection plate please contact your sales rep for more information.

### 96 WELL LOOSE COLUMN RACKS (1ml tubes)

Part Number	Description	Description 2	Guarantee Compatible Collection Plates*	Packaging
30137226	96 Place Rack	Partial plate Cerex 1 cc rack (green)	Greiner	Ea.
253-0006	96 Place Rack	Cerex Only for loose/partial plate work (base plate anodized gray)	Greiner	Ea.
253-0037	96 Place Rack	Cerex Only for loose/partial plate work (base plate anodized gray)	Greiner	Ea.

\*Note: Spacers and rack can also fit other brand of collection plate please contact your sales rep for more information.

## WASTE TRAY AND RACK FOR M10 PROCESSOR

Part Number	Description	Packaging
278-0300	Waste Tray for M10 96	Ea.
279-0013	Waste Tray Rack for M10 48	3 / pk
253-2003	Luer Seal Mat with pull Tab 96 Well	5 / pk

\***Note:** Spacers and rack can also fit other brand of collection plate please contact your sales rep for more information.

## 48 LOOSE TUBE SPACER AND RACKS (1, 3, and 6 ml)

Part Number	Description	Description 2	Packaging
253-0035	ALD3 Luer Seal Rack, 96-place	Facilitates placement of Luer seal on columns	Ea.
279-0014	Collection Rack	For 10x75mm tubes for 48 place system	Ea.
279-0015	Collection Rack	For 12x75mm tubes for 48 Place system	Ea.
279-0016	Collection Rack	For 16x100 tubes for 48mm Place System	Ea.
279-0017	Collection Rack	For 12x32mm Autosampler vials for 48 place systems	Ea.
279-0021	SPE column Rack for 1ml columns	For 48 Place system Flangeless version	Ea.
279-0022	SPE column Rack for 3ml columns	For 48 Place system Flangeless version (if using none flangeless tubes will not fit all 48)	Ea.
279-0023	SPE column Rack for 6ml columns	For 48 Place system Flangeless version (if using none flangeless tubes will not fit all 48)	Ea.
253-0029	Luer Seal Rack	Used to apply Luer seal mat to SPE columns (48)	Ea.

## 96 WELL COLLECTION PLATES

Part Number	Description	Description 2	Packaging
279-0013	96 Well Collection Plate	Greiner 1.2ml Deep Well, Hemispherical	50 / pk
278-0012-SP	96 Well Collection Plate	Greiner 1.2ml Deep Well, Hemispherical	Ea.
278-0060	96 Well Plate Seals	Pre-slit	10 / pk

## VIALS AND CAP

Part Number	Description	Description 2	Packaging
700-2001	Vial Screw Top	12*32mm 9mm	100/PK
700-2002	Vial Screw Top	12*32mm 9mm Conical	100/PK
700-2003	Vial Crimp	12*32mm Conical 11mm	100/PK
700-2004	Vial Crimp	12*32 2ml 11mm	100/PK
700-2005	Vial Screw Top	12*32mm 2ml HC 9MM	100/PK
700-2006	Vial Snap Cap	12*32mm 11mm	100/PK
700-2008	Vial Snap Top	High Rec. 11mm	100/PK

# Simplify your sample preparation.

## REAGENTS

Part Number	Description	Description 2	Packaging
30138399	Reagent Specialized	PFPOH DR 135-100	100g
30142695	Reagent	Silylation MS TFA	100g
30142862	Reagent MTBSTFA	+1% TBDMCS DR102-5	5g
30151125	Reagent	Beta-Gluc. High Efficiency	20ml
700-5007	Reagent DR104	Acylation PFPA	25g
700-5008	Reagent	Acylation HFBA	100g
700-5011	Reagent DR110	Meth. ELUT TMPAH	10ml
700-5014	Reagent DR121	Silylation BSTFA	25g
700-5015	Reagent DR129	Acylation TFAA	100g
700-5018	Reagent Specialized	HFIP	10g
700-5052	Reagent DR2100	Beta-Gluc. 2M	
700-5054	Reagent DR2102	Beta-Gluc 100K Units/ml	25ml
700-5055	Reagent DR2102	Beta-Gluc 100K Units/ml	50ml
700-5057	Reagent	Beta-Gluc 10K Units	10ml
700-5061	Reagent	Beta-Gluc. Recom. E.Coli	20ml
BG100-25ML	Reagent	Beta-Gluc. From Red Abalone	25ml
BG2.5-2MU	Reagent	Beta-Gluc. Lyophilized Power	2mu

+1 626 962 0010

customerservice@tecan.com

**www.tecan.com/lcms**

.....  
**Australia** +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170  
**Italy** +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 595 25 31  
**Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11  
.....

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication; however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

Tecan, Resolvex and Cerex are registered trademarks of Tecan Group Ltd., Männedorf, Switzerland.

© 2019 Tecan Trading AG, Switzerland, all rights reserved.

**www.tecan.com**

