

**Using Consumables Monitoring Disabled Thermo Scientific™ Dionex™ Suppressors  
with  
Thermo Scientific™ Dionex™ ICS-6000 HPIC™ and Integrion™ HPIC™ Systems**

Dear Customer,

In order to manage a global chronic shortage of 1-wire 20k EEPROMS used on our suppressors, we have temporarily removed the consumables monitoring feature from the following suppressors part numbers (P/N).

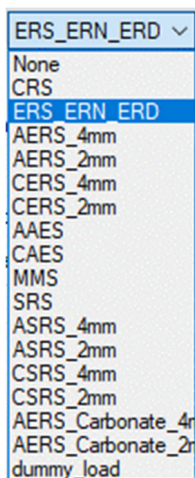
Removal of the consumables monitoring device will have no impact to users not using this feature. Customers utilizing the consumables monitoring feature will not be able to:

- Track the usage of the consumable when moved from system to system
- Run in the Constant Voltage mode

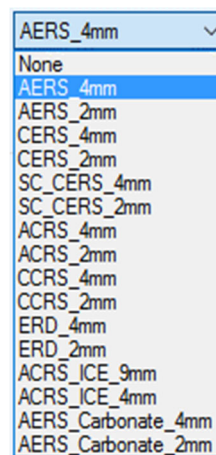
Part number	Suppressor description	P/N of the Consumables Monitoring Disabled (CMD) Suppressors
088668	Thermo Scientific™ Dionex™ CDRS 600 4mm	088668CMD
088670	Thermo Scientific™ Dionex™ CDRS 600 2mm	088670CMD
088666	Thermo Scientific™ Dionex™ ADRS 600 4mm	088666CMD
088667	Thermo Scientific™ Dionex™ ADRS 600 2mm	088667CMD
302663	Thermo Scientific™ Dionex™ CERS 500e 4mm	302663CMD
302664	Thermo Scientific™ Dionex™ CERS 500e 2mm	302664CMD
302661	Thermo Scientific™ Dionex™ AERS 500e 4mm	302661CMD
302662	Thermo Scientific™ Dionex™ AERS 500e 2mm	302662CMD
085087	Thermo Scientific™ Dionex™ ERD 500 4mm	085087CMD
085089	Thermo Scientific™ Dionex™ ERD 500 2mm	085089CMD
083549	Thermo Scientific™ Dionex™ SC-CERS 500 4mm	083549CMD
083547	Thermo Scientific™ Dionex™ SC-CERS 500 2mm	083547CMD

To operate your consumables-monitoring-disabled suppressor with a new instrument method using Dionex ICS-6000 HPIC or Dionex Integrion HPIC systems, please review the menu selections below.

**Expected drop-down menu on the Dionex ICS-6000 HPIC System**



**Expected drop-down menu on the Dionex Integrion HPIC System**



Select the desired suppressor type from the drop-down menu. See the table below for the suggested CMD suppressor for each consumables-monitoring-disabled device.

Suppressor name	CMD Suppressor Part Number	Selection for Dionex ICS-6000 HPIC system	Selection for Dionex Integrion HPIC system
Thermo Scientific Dionex CDRS 600 4mm	088668CMD	CERS_4mm	CERS_4mm
Thermo Scientific Dionex CDRS 600 2mm	088670CMD	CERS_2mm	CERS_2mm
Thermo Scientific Dionex ADRS 600 4mm	088666CMD	AERS_4mm	AERS_4mm
Thermo Scientific Dionex ADRS 600 2mm	088667CMD	AERS_2mm	AERS_2mm
Thermo Scientific Dionex CERS 500e 4mm	302663CMD	CERS_4mm	CERS_4mm
Thermo Scientific Dionex CERS 500e 2mm	302664CMD	CERS_2mm	CERS_2mm
Thermo Scientific Dionex AERS 500e 4mm	302661CMD	AERS_4mm	AERS_4mm
Thermo Scientific Dionex AERS 500e 2mm	302662CMD	AERS_2mm	AERS_2mm
Thermo Scientific Dionex ERD 500 4mm	085087CMD	AERS_4mm	ERD_4mm
Thermo Scientific Dionex ERD 500 2mm	085089CMD	AERS_2mm	ERD_2mm
Thermo Scientific Dionex SC-CERS 500 4mm	083549CMD	CERS_4mm	SC_CERS_4mm
Thermo Scientific Dionex SC-CERS 500 2mm	083547CMD	CERS_2mm	SC_CERS_2mm