

Technical Bulletin

Capillaries for the Thermo Scientific Model 42*i* NO, NO₂, and NO_x Analyzers

Bulletin #112911CAP

Rev 11/2011

The Model 42*i* NO, NO₂, and NO_x Analyzer manuals may contain discrepancies regarding the correct capillary sizes for specific models or configurations. The instrument manuals affected include the Models 42*i*, 42*i*-LS, 42*i*-HL, 42*i*-TL, 42*i*-Y and 42*i*-D.

This technical bulletin contains the most current information regarding correct capillary sizes. When performing routine maintenance, please ensure that the correct size capillaries are installed per the table below.

Please note that the manual discrepancies do not impact any instruments at the time of shipment. All 42*i* series instruments were built and shipped with the correct capillaries as listed below.

The next revisions to the instrument manuals will reflect the correct capillary size.

Model	Converter	Pump Type	Sample		Ozone		Bypass		Oxygen ²	
			P/N	Size ¹	P/N	Size ¹	P/N	Size ¹	P/N	Size ¹
42 <i>i</i>	Molybdenum	Internal	4127	15	4119	8	NA	NA	NA	NA
42 <i>i</i>	Stainless Steel ³	Internal	4121	10	4119	8	NA	NA	NA	NA
42 <i>i</i> -LS	Moly or SS	Internal	4119	8	4121	10	NA	NA	4119	8
42 <i>i</i> -HL	Moly or SS	External	4113	5	4121	10	4126	20	4119	8
42 <i>i</i> -TL	Molybdenum	External	4126	20	4119	8	NA	NA	NA	NA
42 <i>i</i> -Y	Molybdenum	External	4126	20	4119	8	4110	22	NA	NA
42 <i>i</i> -D	Moly or SS	Internal	4127	15	4119	8	NA	NA	NA	NA

¹ Size in milli-inches (mil)

² Only when option is installed

³ When Stainless Steel Converter (SS) is installed, sample flow is 0.200 LPM and flow alarms should be set as follows: 0.150 LPM (Min) and 0.300 LPM (Max)

Capillary Reference Guide

Part Number	Capillary Size	Color
4110	22 milli-inches	Yellow / Black
4113	5 milli-inches	Yellow
4119	8 milli-inches	Red
4121	10 milli-inches	Black
4126	20 milli-inches	Blue
4127	15 milli-inches	Purple

If you have any questions or require technical assistance, please contact your local distributor or our Technical Support team via telephone at (508) 520-0430 or by e-mail at AQIService@thermofisher.com.