

Easy-fit ISIS 3 Sample Loops recommended sample introduction timing

Using our [Easy-fit fixed-volume sample loops](#), you can benefit from these default timing recommendations that we have determined so you can get your method set up and start your analysis quicker.

These are our recommended timings and pump speeds for each sample loop in the range. These are based upon our measurements with an Agilent 7900 ICP-MS configured with ISIS 3 7-port valve and SPS 4 autosampler fitted with standard tubing:

- 1.02mm ID white/white peristaltic pump tubing for carrier [5005-0020](#)
- 0.25mm ID blue/orange peristaltic pump tubing for Internal Standard (ISTD)/Tune [5005-0021](#)
- 1mm ID sample probe [G8410-80103](#)

Part Number	Loop Tubing ID (mm)	Loop Volume (mL)	Typical Acquisition Time (sec)	Sample Load (sec) @ Speed 60% ISIS Uptake Pump	Stabilization Time (sec)	Loop Probe Wash (sec) @ Speed 80%	Loop Wash (sec) @ Speed 80%
5005-0030	1	0.25	15	7	20	4	7
5005-0031	1	0.50	25	8	20	4	8
5005-0032	1	0.75	50	9	20	4	9
5005-0033	1	1.00	80	10	20	4	10
5005-0034	1	1.25	115	11	20	4	11
5005-0035	1	1.50	145	12	20	4	12
5005-0036	1	2.00	210	14	20	4	14
5005-0037	1	2.50	375	17	20	4	17
5005-0038	2.18	1.00	80	8	20	4	8
5005-0039	2.18	1.25	105	9	20	4	9
5005-0040	2.18	1.50	130	10	20	4	10
5005-0041	2.18	2.00	260	11	20	4	11
5005-0042	2.18	2.50	365	12	20	4	12
5005-0043	2.18	3.00	410	14	20	4	14
5005-0044	2.18	4.00	645	17	20	4	17

Recommended timings are based on default settings; optimization for your application and system may be required. Please refer to the ISIS 3 user guide for further information.

www.agilent.com/chem/easyfitloops

DE74080043

This information is subject to change without notice.