

Sample name:	Ciclotizolam	Operator :	Jianmei Liu
Data file:	37999-0668803.dx	Instrument:	ISO GCMS2
Acq. method:	GC-MSD General Method 3.amx	Inj. volume:	1.000
Acquired on:	2023-05-12 11:21:15-04:00	Column:	
Method Description:			

