

Sample name: Flubromazepam
Data file: 15157-0570041.dx
Acq. method: GC-MSD General Method 2.amx
Acquired on: 2020-01-30 19:38:56-05:00
Method Description:

Operator : Samantha Goodwin
Instrument: ISO GCMS2
Inj. volume: 1.000
Column:

