

Sample name: 2-fluoro-2-oxo PCE (hydrochloride)

Data file: 36722-0644444.dx

Operator : Jianmei Liu

Acq. method: GC-MSD General Method 1.amx

Instrument: ISO GCMS2

Acquired on: 2022-09-26 15:01:54-04:00

Inj. volume: 1.000

Method Description:

Column:

