

Sample name: N-cyclohexyl Butylone (hydrochloride)

Data file: 36884-0646340.dx

Operator : Jianmei Liu

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

Instrument: ISO GCMS2

Acquired on: 2022-09-26 15:56:01-04:00

Inj. volume: 1.000

Method Description:

Column:

