

Sample name: 9(R)-Hexahydrocannabinol

Data file: 27500-0646274.dx

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

Acquired on: 2022-06-09 12:02:35-04:00

Method Description:

Operator : Jianmei Liu

Instrument: ISO GCMS2

Inj. volume: 1.000

Column:

