

Sample name: Cannabidiol diacetate
Data file: 36424-0646266.dx
Acq. method: GC-MSD General Method 1.amx
Acquired on: 2023-02-25 09:35:31-05:00
Method Description:

Operator : Monica Bugeja
Instrument: ISO GCMS2
Inj. volume: 1.000
Column:

