

File :C:\ISO\19APR2018-NON ISO\10979-0528272.D
Operator : JianmeiL
Acquired : 20 Apr 2018 2:30 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 3,4-MDMA methylene homolog (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 19

