

File :Y:\11047-0455466.D
Operator :TomB
Acquired : 27 May 2015 10:47 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: JWH 018 N-(5-bromopentyl) analog
Misc Info :
Vial Number: 19

