

File : C:\MassHunter\Data\2024\2024-12-06 non ISO\38176-0689195.D
Operator : Jianmeil/ Camille W
Acquired : 07 Dec 2024 02:54 using AcqMethod GC MS General method 1.M
Instrument : ISO GC MS3
Sample Name: Pyrazolam-d7
Misc Info : 50C hold 1 min 30C/min to 300C
Vial Number: 23

