

Sample name: 1-(2-Chloro-3-fluorophenyl)piperazine (hydrochloride)
Data file: 29116-0572379.dx **Operator :** Samantha Goodwin
Acq. method: GC-MSD General Method 2.amx **Instrument:** ISO GCMS2
Acquired on: 2020-01-30 17:19:45-05:00 **Inj. volume:** 1.000
Method Description: **Column:**

