

³Thermo Fisher Scientific, Winnipeg, MB, Canada

