

dataTaker is pleased to announce the use of 3G is available in the DT8xM range of product for the US market.

What this means for US customers

The 3G modem supplied within the DT8xM range is FCC and PTCRB tested for use on US networks but is not AT&T approved.

Therefore the DT8xM 3G models will be offered for sale to be used on other US networks such as T-Mobile.

An upgrade of existing DT8xM 2G models with the 3G modem will be made available and details will be realised to distributors in due course.

For further details contact your local Thermo Fisher Scientific sales professional.





05-06-2014