

## Monolith Explorer Kit – Mission to Mars (Cat# MO-X001)

Single labeling reaction with RED-NHS and RED-tris-NTA 2nd Generation dyes.

The goal of this mission is to get your assay development going before you commit to a full kit. We've packed everything you'll need to perform one independent labeling reaction for proteins and peptides with RED-NHS and RED-tris-NTA 2nd Generation dyes. Includes dyes, dye removal reagents, and 60 premium capillaries.

Included in this kit	Quantity	Storage
Monolith NT.115 premium capillaries	60	Dry and dark at RT
Dye RED-NHS 2nd Generation (10 µg)	1	-20 °C
Labeling buffer NHS	1	4 °C
A-Column	1	4 °C
B-Column	1	4 °C
Column adapter	1	RT
Product insert (MO-L011)	1	RT
Dye RED-tris-NTA 2nd Generation (125 pmol)	1	-20 °C
5x PBS-T (2 mL)	1	-20 °C
His6 control peptide (1200 pmol)	1	-20 °C
Product insert (MO-L018)	1	RT

### Capillary handling and safety information

Keep vial closed to ensure capillaries are clean and dust-free. Hold individual capillaries at one end, not in the center. Use gloves. Dispose of capillaries according to applicable regulations and according to the sample contained in them.



**WARNING** Broken glass can cut skin. Handle capillaries carefully to avoid breaking. Do not touch broken capillaries.

### Packaging disposal

All packaging, including the cardboard insert with foil, can be recycled with paper products.

### Contact



Need help? Visit the NanoTemper Explorer Community at [nanotempertech.com/explore](https://nanotempertech.com/explore)

