

MB-TCO

A commonly used redox indicator in nucleic acid research. MB-TCO (trans-cyclooctene) may be used to label tetrazine tagged molecules via a copper-free reaction.



Product attributes

Product Description

Methylene Blue (MB) is a commonly used redox indicator in nucleic acid research. It is also studied for its use in medical applications as well as being used as a general biological stain. Reactive formats of MB can be conjugated to biomolecules. The conjugate will have a blue color and be able to complex with nucleic acids.

MB-TCO (trans-cyclooctene) is a reactive form of MB that may be used to label tetrazine tagged molecules via a copper-free reaction. We also offer [MB Acid](#) and a selection of other chemically reactive formats for use in labeling biomolecules such as proteins and nucleic acids.

- MW: ~650
- Store at -20 °C and protected from light

See the table below for our full list of methylene blue derivatives and formats.

Methylene Blue Derivatives

| Product | Size | Catalog No. | Features |
|---------------------------------------|------|-----------------------|--|
| MB Acid | 5 mg | 40076 | Free acid form |
| MB Succinimidyl Ester | 5 mg | 40075 | Amine-reactive chemistry for labeling proteins |
| MB-Maleimide | 1 mg | 40118 | Thiol-reactive chemistry for labeling proteins |
| MB-DBCO | 1 mg | 40114 | Allows bioorthogonal conjugation to label azide containing molecules |
| MB-Methyltetrazine | 1 mg | 40115 | Allows labeling of TCO tagged molecules |
| MB-TCO | 1 mg | 40116 | Allows labeling of tetrazine tagged molecules |
| MB-Azide | 1 mg | 40117 | Allows labeling alkyne, BCN, or phosphine-containing molecules. |

See our other [reactive DNA/RNA binding dyes](#).

This datasheet was generated on January 21, 2026 at 06:24:58 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mb-tco/>