

## MB Succinimidyl Ester

The amine reactive succinimidyl ester of MB (methylene blue) can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. The conjugate will have a blue color and be able to complex with nucleic acids.



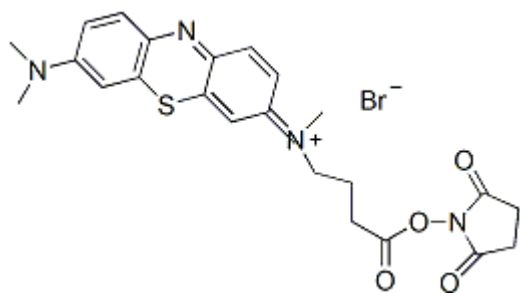
### Product attributes

## 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.

The amine reactive MB Succinimidyl Ester can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. We also offer [MB Acid](#) and a selection of other chemically reactive formats for use in labeling biomolecules such as proteins and nucleic acids.

- $C_{24}H_{26}BrN_4O_4S$
- Blue solid soluble in DMF or DMSO
- Store at  $-20^{\circ}C$ , desiccated
- MW: 533



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

## Methylene Blue Derivatives

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

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