

www.biotium.com

Revised: January 8, 2023

Product Information

Methylene Blue (MB) Derivatives

Product List

Cat. No.	Product	Unit Size	MW
40075	MB Succinimidyl Ester	5 mg	533
40076	MB Acid	5 mg	392
40114	MB-DBCO	1 mg	~790
40115	MB-Methyltetrazine	1 mg	~620
40116	MB-TCO	1 mg	~650
40117	MB-Azide	1 mg	~520
40118	MB-Maleimide	1 mg	~560

Storage and Handling

Store MB Acid at 4°C and protected from light; store all other MB derivatives at -20°C and protected from light. Store MB Succinimidyl Ester and MB-Maleimide solids desiccated. Product is stable for at least 12 months from date of receipt when stored as recommended.

Prepare the stock solution in DMSO. The MB products are not soluble in other solvents. Solutions can be aliquoted and stored with desiccant and protected from light at -20°C, for up to 6 months.

Spectral Properties

Methylene Blue shows a primary absorption peak at 665 nm with a shoulder around 610 nm, and a small peak at around 293 nm [1].



Figure 1. Left: MB Succinimidyl Ester, Right: MB Acid, (others proprietary).

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. We offer MB Free Acid and with a selection of six different chemically reactive groups for use in labeling biomolecules such as proteins and nucleic acids. For more conventional labeling chemistry, we offer MB Succinimidyl Ester, MB-Maleimide, and MB Acid (Figure 1). MB Succinimidyl Ester is reactive with primary amines and can be used to label proteins or other molecules with primary amine groups. MB Maleimide can be used to react with free thiol groups. The resulting conjugates will have a blue color and be able to complex with nucleic acids. MB Acid can be used as a non-reactive control dye for conjugates.

We offer four MB conjugates for bioorthogonal labeling reactions (precise structures proprietary). MB-dibenzocyclooctyne (DBCO) allows copper-free bioorthogonal conjugation to spontaneously label molecules containing azide groups. MB-methyltetrazine and MB-TCO (trans-cyclooctene) react with the corresponding TCO, tetrazine and methyltetrazine via a copper-free reaction. MB-Azide allows a copper-catalyzed bioorthogonal conjugation reaction with alkyne, copper-free reaction with BCN, or Staudinger ligation with phosphine. We do not provide a validated protocol for these products and ask that customers use the related literature to find appropriate guidance for their own application.

Biotium also offers reactive formats of fluorescent CF® Dyes for labeling proteins, nucleic acids, or other biomolecules (see Related Products).

References

1. J Clin Med, 9, 11, (2020).

Cat. No.	Product
40020-1	5-Aminoallyl-dUTP (AA-dUTP), Lyophilized Powder
40020	5-Aminoallyl-dUTP (AA-dUTP), 10 mM solution
40021	5-Aminoallyl-UTP, Sodium Salt, 10 mM Solution
40021-1	5-Aminoallyl-UTP, Sodium Salt, Lyophilized Powder
90117-90119	Cyanine NS Dye Succinimidyl Ester
9208096000	CF® Dye Azide
9218796007	CF® Dye Picolyl Azide
92086-92090	CF® Dye Alkyne
9210397502	CF® Dye SE/TFP
9202096079	CF® Dye Maleimides
96028-96039	CF® Dye Methyltetrazine
96040-96051	CF® Dye TCO
92167	Biotin Azide
92168	Biotin Alkyne
92169	Biotin BCN
92186	Biotin Picolyl Azide
92100-92101	Cyanine Dye Alkyne
96023	MTS-BCN
96038	Biotin Methyltetrazine
96050	Biotin TCO

Please visit our website at www.biotium.com for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF® Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

CF® Dye technology is covered by U.S. and international patents

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.