

**Buccutite™ MTA, maleimide [MTAM]**Catalog number: 5358  
Unit size: 2 umoles**Product Details**

---

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

**Chemical Properties**

---

Appearance	Solid
Molecular Weight	514.54
Soluble In	DMSO

**Applications**

---

Buccutite™ crosslinking technology provides the most convenient and effective crosslinking method to link two biomolecules with a high conjugation yield. The method uses one pair of crosslinkers: Buccutite™ MTA and Buccutite™ FOL. MTA is added to one molecule, while FOL is added to another molecule. The cross-linking reaction is initiated by mixing Molecule-1-Buccutite™ MTA and Molecule-2-Buccutite™ FOL under neutral conditions. Many of our customer have requested us to offer the stand-alone Buccutite™ MTA and Buccutite™ FOL reagents to expand the application of Buccutite™ crosslinking technology. Buccutite™ MTA maleimide (MTAM) can be used the same way as the widely used SMCC for crosslinking proteins. One end of the MTAM reacts (via maleimide) with thiols (-SH) of cysteine found in the reduced antibodies (by TCEP or DTT). SMCC crosslinking requires high concentration of proteins. In addition, SMCC-modified protein is extremely unstable and often self-reactive since proteins often contain both amine and thiol groups that cause significant amount of homo-crosslinking. Buccutite™ crosslinking reaction occurs under extremely mild and neutral conditions without any catalyst required. It is robust and efficient.