

Product Information

CPMTS

Catalog #: 91011

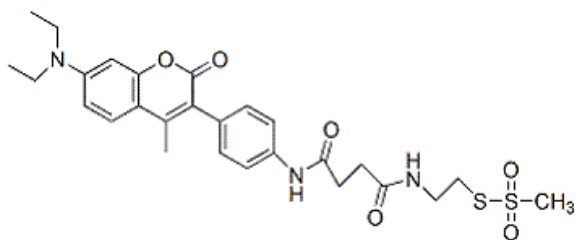
Unit Size: 5 mg

Molecular Formula: C₂₇H₃₃N₃O₆S₂

Molecular Weight: ~560

Color & Form: Yellow solids

Structure:



Storage and Handling

Store CPMTS at -20°C and protected from light. Stock solution can be prepared in DMSO or DMF and can be stored at ≤ -20°C for at least 12 months.

Product Description

Biotium developed CPMTS as an alternative to the widely used thiol-reactive coumarin dye CPM (#91010). CPMTS and CPM have nearly identical absorption and emission spectra. However, CPMTS contains a MTS group, which could make the dye more advantageous than CPM for sulfhydryl labeling because of the high selectivity and reactivity of MTS reagents toward thiol groups.

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