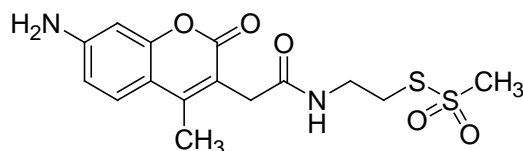


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: MTS-AMCA**CATALOG #** 91033**MOLECULAR INFORMATION:** C₁₅H₁₈N₂O₆S₂
MWt: 370.45**PROPERTIES:**

Color & Form Of-white solid
Purity > 98% by TLC
Solubility DMF or DMSO
Absorption/Emission 353/442 nm (MeOH)
Extinction Coefficient

STORAGE AND HANDLING: Store at -20 °C and protect from light

APPLICATION: AMCA (7-amino-4-methylcoumarin-3-acetic acid) is one of the most widely used UV-excitable, blue fluorescent dye. MTS-AMCA readily labels a thiol with AMCA.

TOXICITY: Not established. Not listed by NTP, IARC or OSHA.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--