

MTS-AMCA

MTS-AMCA (7-amino-4-methylcoumarin-3-acetic acid methanethiosulfonate) readily labels a thiol with AMCA, a widely used UV-excitable, blue fluorescent dye.



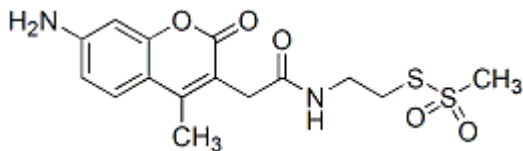
Product attributes

Excitation/Emission	353/442 nm
---------------------	------------

Product Description

MTS-AMCA (7-amino-4-methylcoumarin-3-acetic acid methanethiosulfonate) readily labels a thiol with AMCA, a widely used UV-excitable, blue fluorescent dye.

- $\lambda_{Ex}/\lambda_{Em}$ = 353/442 nm
- Off-white solid soluble in DMF or DMSO
- Store at -20 °C and protect from light
- $C_{15}H_{18}N_2O_5S_2$
- MW: 370.45



This datasheet was generated on June 22, 2026 at 01:12:36 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mts-amca/>