



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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.



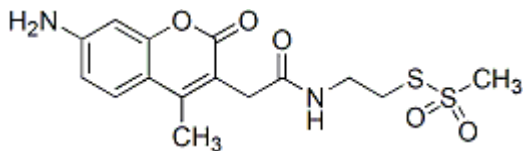
Product attributes

| | |
|---------------------|------------|
| 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 December 17, 2025 at 09:39:25 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mts-amca/>