

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

## PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 6-((7-amino-4-methylcoumarin-3-acetyl) amino)hexanoic acid,

succinimidyl ester (AMCA-X SE)

**CATALOG #** 90077

**MOLECULAR**  $C_{22}H_{25}N_3O_7$  **INFORMATION:** Mwt: 443.46

$$\begin{array}{c|c} H_2N & O & O & O \\ \hline & O & O & O \\ \hline & NH(CH_2)_5C-O-N \\ \hline & O & O \\ \hline \end{array}$$

**PROPERTIES:** 

**Color & Form** Off-white solid

**Solubility** Soluble in DMSO or DMF **Absorption/Emission** 353/442nm (MeOH)

Extinction 19,000

**STORAGE AND** Store desiccated at -20 °C. Protect from light, especially in solution. **HANDLING:** 

**APPLICATION:** AMCA-X SE is an amine-reactive blue fluorescent dye widely used for labeling

proteins or other biomolecules that contain a primary or secondary aliphatic amine. A stock solution of the dye can be made in anhydrous DMF or DMSO. The coupling reaction is usually carried out in a pH ~8.5 buffer (for example, 0.1-0.2 M sodium bicarbonate buffer). Alternatively, the coupling reaction may also be carried out in an organic solvent for amino compounds that are insoluble in aqueous solution.

Ref. 1) J Immunol **164**, 5236(2000); 2) J Biol Chem **276**, 10366(2001).

**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:** 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.