

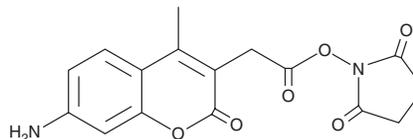
# PRODUCT INFORMATION



## AMCA N-succinimidyl ester

Item No. 34668

**CAS Registry No.:** 113721-87-2  
**Formal Name:** 7-amino-4-methyl-2-oxo-2H-1-benzopyran-3-acetic acid, 2,5-dioxo-1-pyrrolidinyl ester  
**Synonyms:** 7-Amino-4-methylcoumarin-3-acetic Acid N-succinimidyl ester, 7-Amino-4-methyl-3-coumarinylacetic Acid N-succinimidyl ester  
**MF:** C<sub>16</sub>H<sub>14</sub>N<sub>2</sub>O<sub>6</sub>  
**FW:** 330.3  
**Purity:** ≥90%  
**Abs./Em. Max:** UV/400-460 nm  
**UV/Vis.:** λ<sub>max</sub>: 236, 359 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

AMCA N-succinimidyl ester is supplied as a solid. A stock solution may be made by dissolving the AMCA N-succinimidyl ester in the solvent of choice, which should be purged with an inert gas. AMCA N-succinimidyl ester is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AMCA N-succinimidyl ester in these solvents is approximately 30 mg/ml.

AMCA N-succinimidyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AMCA N-succinimidyl ester should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AMCA N-succinimidyl ester has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

AMCA N-succinimidyl ester is an amine-reactive fluorescent probe and ester form of AMCA (Item No. 29184).<sup>1</sup> It contains an N-hydroxysuccinimide ester that reacts with lysine residues to form photostable amide links. Upon activation with UV light, AMCA displays emission maxima of 400-460 nm. AMCA N-succinimidyl ester has commonly been used in multiplex immunophenotyping applications.<sup>2-4</sup>

### References

1. Villalobos-Molina, R., Gallardo-Ortíz, I.A., López-Guerrero, J.J., *et al.* Evidence that NAN-190-induced hypotension involves vascular α<sub>1</sub>-adrenoceptor antagonism in the rat. *Eur. J. Pharmacol.* **455(1)**, 59-64 (2002).
2. Ferri, G.-L., Gaudio, R.M., Castello, I.F., *et al.* Quadruple immunofluorescence: A direct visualization method. *J. Histochem. Cytochem.* **45(2)**, 155-158 (1997).
3. Nederlof, P.M., Robinson, D., Abuknesha, R., *et al.* Three-color fluorescence in situ hybridization for the simultaneous detection of multiple nucleic acid sequences. *Cytometry* **10(1)**, 20-27 (1989).
4. Gothot, A., Grosdent, J.-C., and Paulus, J.-M. A strategy for multiple immunophenotyping by image cytometry: Model studies using latex microbeads labeled with seven streptavidin-bound fluorochromes. *Cytometry* **24(3)**, 214-225 (1996).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/22/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM