520 Mercury Drive, Sunnyvale, CA 94085 Tel: 408-733-1055; Fax: 408-733-1304

www.aatbio.com; E-Mail: info@aatbio.com

Product Information Sheet

Ordering Information

Product Number: 551

Product Name: 7-Hydroxycoumarin-3-carboxylic acid, succinimidyl ester

Unit Size: 50 mg

Storage Conditions: <-15 °C and kept from light and moisture

Expiration Date: 12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 303.23

Chemical Structure:

Appearance: Off-white solid

Soluble in: DMSO
Excitation Wavelength: 419 nm
Emission Wavelength: 447 nm

Application Notes

7-Hydroxycoumarin-3-carboxylic acid, SE is the amine-reactive succinimidyl ester of 7-Hydroxycoumarin-3-carboxylic acid. It is a blue fluorescent dyes for labeling proteins and nucleic acids. This coumarin is also increasingly used to label peptides, nucleotides and carbohydrates.