

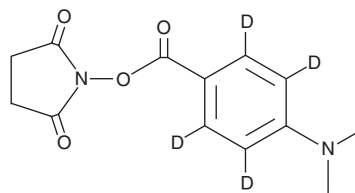
# PRODUCT INFORMATION



## DMABA-d<sub>4</sub> NHS ester

Item No. 11217

**CAS Registry No.:** 1175002-03-5  
**Formal Name:** 4-(dimethylamino)-benzoic-2,3,5,6-d<sub>4</sub> acid, 2,5-dioxo-1-pyrrolidinyl ester  
**Synonyms:** 4-(dimethylamino) Benzoic Acid-d<sub>4</sub>, DMABA N-hydroxysuccinimide ester-d<sub>4</sub>, N-Succinimidyl 4-(dimethylamino)benzoate-d<sub>4</sub>  
**MF:** C<sub>13</sub>H<sub>10</sub>D<sub>4</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 266.3  
**Chemical Purity:** ≥98% (DMABA NHS ester\_)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>4</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 234, 322 nm  
**Supplied as:** A crystalline 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

DMABA-d<sub>4</sub> NHS ester is intended for use in combination with DMABA NHS ester (Item No. 11216) as a derivatizing reagent for Phosphatidylethanolamine (PE) lipids in order to facilitate MS characterization and to quantify relative changes in their abundance.

DMABA-d<sub>4</sub> NHS ester is supplied as a crystalline solid. A stock solution may be made by dissolving the DMABA-d<sub>4</sub> NHS ester in the solvent of choice, which should be purged with an inert gas. DMABA-d<sub>4</sub> NHS ester is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of DMABA-d<sub>4</sub> NHS ester in these solvents is approximately 20 mg/ml.

### Description

DMABA-d<sub>4</sub> and DMABA NHS esters have been shown to facilitate the characterization of PE lipids via electrospray mass spectrometry by providing a common precursor ion scan for diacyl, ether, and plasmalogen PE lipids.<sup>1</sup> Furthermore, application of split-and-pool methodology allows for relative changes in endogenous PE lipid levels between control and treated samples to be measured.<sup>2</sup> PE lipids are important components of cell membranes and biochemical pathways of fatty acid synthesis that contain abundant polyunsaturated fatty acyl groups. Oxidation of these phospholipids may be linked to various human diseases. DMABA NHS ester is a reagent that reacts with the primary amine group of PE lipids.<sup>1,2</sup> This facilitates the use of electrospray tandem mass spectrometry for the detection of diacyl, ether, and plasmalogen PE lipids that cannot be readily observed otherwise.<sup>1,2</sup>

### References

1. Zemski Berry, K.A., Turner, W.W., VanNieuwenhze, M.S., et al. Stable isotope labeled 4-(dimethylamino) benzoic acid derivatives of glycerophosphoethanolamine lipids. *Anal. Chem.* **81(16)**, 6633-6640 (2009).
2. Zemski Berry, K.A., Turner, W.W., VanNieuwenhze, M.S., et al. Characterization of oxidized phosphatidylethanolamine derived from RAW 264.7 cells using 4-(dimethylamino)benzoic acid derivatives. *Eur. J. Mass Spectrom* **16(3)**, 463-470 (2010).

#### 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, 02/22/2024

#### 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