

# PRODUCT INFORMATION



## L-ANAP methyl ester (hydrochloride)

Item No. 15437

**Formal Name:** 3-[(6-acetyl-2-naphthalenyl)amino]-L-alanine, methyl ester, monohydrochloride

**MF:** C<sub>16</sub>H<sub>18</sub>N<sub>2</sub>O<sub>3</sub> • HCl

**FW:** 322.8

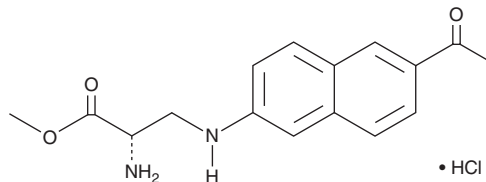
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 248, 256, 354 nm

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥2 years



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

### Laboratory Procedures

L-ANAP methyl ester (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the L-ANAP methyl ester (hydrochloride) in the solvent of choice, which should be purged with an inert gas. L-ANAP methyl ester (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of L-ANAP methyl ester (hydrochloride) in these solvents is approximately 0.5, 16, and 14 mg/ml, respectively.

### Description

L-ANAP methyl ester is an esterified form of the fluorescent unnatural amino acid L-ANAP (Item No. 15436).<sup>1</sup> It has been used in the synthesis of L-ANAP. L-ANAP methyl ester has also been used to fluorescently label proteins *in vitro* through genetic encoding.<sup>2</sup>

### References

1. Xiang, Z. and Wang, L. Enantiospecific synthesis of genetically encodable fluorescent unnatural amino acid L-3-(6-acetylnaphthalen-2-ylamino)-2-aminopropanoic acid. *J. Org. Chem.* **76(15)**, 6367-6371 (2011).
2. Berlin, J.R., Lopez, W., Jain, M.R., *et al.* Use of the methyl ester of a fluorescent unnatural amino acid to facilitate site-specific incorporation of fluorescent probes in proteins. *Biophys. J.* **108(2)**, 624A-625A (2015).

#### 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, 04/07/2022

#### 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