PRODUCT INFORMATION



L-ANAP (trifluoroacetate salt)

Item No. 15436

CAS Registry No.:	2287346-69-2	0
Formal Name:	3-[(6-acetyl-2-naphthalenyl)amino]-	a a Ŭ
	L-alanine, monotrifluoroacetate salt	
MF:	$C_{15}H_{16}N_2O_3 \bullet CF_3COOH$	0
FW:	386.3	
Purity:	≥95%	
UV/Vis.:	λ _{max} : 227, 247, 350 nm	NH ₂ H
Supplied as:	A crystalline solid	• CE ₂ COOH
Storage:	-20°C	01300011
Stability:	≥4 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

Laboratory Procedures

L-ANAP (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the L-ANAP (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. L-ANAP (trifluoroacetate salt) is soluble in ethanol, DMSO and dimethyl formamide (DMF). The solubility of L-ANAP (trifluoroacetate salt) in these solvents is approximately 0.5 mg/ml in ethanol and approximately 10 mg/mg in DMSO and DMF.

Description

L-ANAP is an unnatural amino acid with intrinsic fluorescence that can be genetically encoded into proteins.¹ This technology allows incorporation of L-ANAP into virtually any site on a protein, providing a unique method for imaging biological processes in vivo.

Reference

1. Lee, H.S., Guo, J., Lemke, E.A., et al. Genetic incorporation of a small, environmentally sensitive, fluorescent probe into proteins in Saccharomyces cerevisiae. J. Am. Chem. Soc. 131(36), 12921-12923 (2009).

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.

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, 12/05/2022

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