PRODUCT INFORMATION



NAB 2

Item No. 26145

CAS Registry No.:	1504588-00-4		0	
Formal Name:	N-[(2-chlorophenyl)methyl]-	\rightarrow \wedge		
	1-(2,5-dimethylphenyl)-1H-		N'	
	benzimidazole-5-carboxamide		ή (I	
MF:	$C_{23}H_{20}CIN_3O$			N'
FW:	389.9			
Purity:	≥98%			
UV/Vis.:	λ _{max} : 220 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

NAB 2 is supplied as a crystalline solid. A stock solution may be made by dissolving the NAB 2 in the solvent of choice, which should be purged with an inert gas. NAB 2 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Description

NAB 2 is an inhibitor of α -synuclein-induced toxicity.¹ It prevents α -synuclein-induced defects in Rsp5-dependent endocytosis and vacuolar delivery of Mup1, Golgi-to-vacuole trafficking of Sna3, and degradation of Bap2 in wild-type yeast cells. NAB 2 also prevents formation of α -synuclein foci in wild-type yeast cells, an effect that is prevented by the *rsp5*^{G747E} point mutation.

Reference

1. Tardiff, D.F., Jui, N.T., Khurana, V., et al. Yeast reveal a "druggable" Rsp5/Nedd4 network that ameliorates a-synuclein toxicity in neurons. Science 342(6161), 979-983 (2013).

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, 10/26/2022

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