PRODUCT INFORMATION



U-93631

Item No. 26417

CAS Registry No.:	152273-12-6	
Formal Name:	4,5-dihydro-4,4-dimethyl-imidazo[1,5-a]	
	quinoxaline-3-carboxylic acid,	
	1,1-dimethylethyl ester	
MF:	$C_{17}H_{21}N_{3}O_{2}$	N
FW:	299.4	Ń I
Purity:	≥98%	
UV/Vis.:	λ _{max} : 222, 325 nm	
Supplied as:	A crystalline solid	
Storage:	-20°C	0
Stability:	≥4 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

Laboratory Procedures

U-93631 is supplied as a crystalline solid. A stock solution may be made by dissolving the U-93631 in the solvent of choice, which should be purged with an inert gas. U-93631 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of U-93631 in ethanol and DMSO is approximately 20 mg/ml and approximately 25 mg/ml in DMF.

Description

U-93631 is a $GABA_A$ receptor ligand.¹ It accelerates decay of GABA-induced currents in HEK293 cells expressing rat $\alpha 1\beta 2$, $\beta 2\gamma 2$, and $\alpha 1\gamma 2$ subunit-containing GABA_A receptors when used at a concentration of 5 μ M. U-93631 inhibits radioligand binding to the picrotoxin site on α 1 β 2 γ 2 subunit-containing GABA_A receptors.²

References

- 1. Dillon, G.H., Im, H.K., Hamilton, B.J., et al. U-93631 causes rapid decay of y-aminobutyric acid-induced chloride currents in recombinant rat γ -aminobutyric acid type A receptors. Mol. Pharmacol. 44(4), 860-865 (1993).
- 2. Dillon, G.H., Im, W.B., Pregenzer, J.F., et al. [4-Dimethyl-3-t-butylcarboxyl-4,5-dihydro (1,5-a) quinoxaline] is a novel ligand to the picrotoxin site on GABA_A receptors, and decreases single-channel open probability. J. Pharmacol. Exp. Ther. 272(2), 597-603 (1995).

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, 12/20/2020

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