PRODUCT INFORMATION



AM3102

Item No. 13452

CAS Registry No.: Formal Name:	213182-22-0 N-[(1R)-2-hydroxy-1-methylethyl- 97-octadecenamide	ц
Synonym:	KDS-5104	Ï
MF:	$C_{21}H_{41}NO_2$	
FW:	339.6	
Purity:	≥98%	
Supplied as:	A crystalline solid	\vee \vee \vee \vee
Storage:	-20°C	
Stability:	≥4 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

Laboratory Procedures

AM3102 is supplied as a crystalline solid. A stock solution may be made by dissolving the AM3102 in the solvent of choice, which should be purged with an inert gas. AM3102 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AM3102 in these solvents is approximately 5, 20, and 25 mg/ml, respectively.

Description

AM3102 is an OEA analog that stimulates PPAR α transcriptional activity with an EC₅₀ value of 100 nM and prolongs feeding latency in rodents with an ED₅₀ value of 2.4 mg/kg.¹ It is resistant to enzymatic hydrolysis and is as potent as OEA in enhancing transcriptional activity of PPAR α and reducing food intake in free-feeding rats.^{1,2} AM3102 demonstrates weak affinity for the central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors with K_i values of 33 and 26 μ M, respectively.³

References

- 1. Astarita, G., Di Giacomo, B., Gaetani, S., et al. Pharmacological characterization of hydrolysis-resistant analogs of oleoylethanolamide with potent anorexiant properties. J. Pharmacol. Exp. Ther. 318(2), 563-570 (2006).
- 2. Wang, J. and Ueda, N. Role of the endocannabinoid system in metabolic control. Curr. Opin. Nephrol. Hypertens. 17(1), 1-10 (2008).
- 3. Lin, S., Khanolkar, A.D., Fan, P., et al. Novel analogues of arachidonylethanolamide (anandamide): Affinities for the CB1 and CB2 cannabinoid receptors and metabolic stability. J. Med. Chem. 41(27), 5353-5361 (1998).

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

uyer 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, 03/21/2024

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