

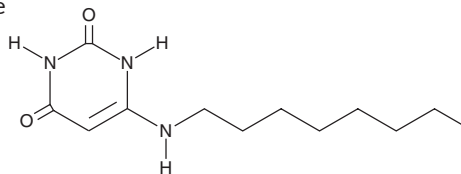
PRODUCT INFORMATION



6-OAU

Item No. 17687

CAS Registry No.: 83797-69-7
Formal Name: 6-(octylamino)-2,4(1H,3H)-pyrimidinedione
MF: C₁₂H₂₁N₃O₂
FW: 239.3
Purity: ≥98%
UV/Vis.: λ_{max}: 265 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

6-OAU is supplied as a crystalline solid. A stock solution may be made by dissolving the 6-OAU in the solvent of choice, which should be purged with an inert gas. 6-OAU is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 6-OAU in these solvents is approximately 0.25 mg/ml.

Description

GPR84 is a Gα_i protein-coupled receptor that is activated by medium-chain free fatty acids, but only at micromolar concentrations.¹ GPR84 is expressed in leukocytes and its expression is further induced by LPS.^{1,2} 6-OAU is a synthetic agonist for GPR84 (EC₅₀ = 512 nM).² It has insignificant actions at a panel of 69 other receptors, channels, or transporters.² 6-OAU promotes chemotaxis in human neutrophils (EC₅₀ = 318 nM) and amplifies TNF-α production by LPS-treated U937 macrophages.² In rats, 6-OAU (1 mg/ml) drives the recruitment of neutrophils and macrophages when injected into a subcutaneous air pouch and raises serum levels of CXCL1 (growth-regulated α protein) when delivered intravenously.²

References

1. Davenport, A.P., Alexander, S.P.H., Sharman, J.L., *et al.* International Union of Basic and Clinical Pharmacology. LXXXVIII. G protein-coupled receptor list: Recommendations for new pairings with cognate ligands. *Pharmacol. Rev.* **65**(3), 967-986 (2013).
2. Suzuki, M., Takaishi, S., Nagasaki, M., *et al.* Medium-chain fatty acid-sensing receptor, GPR84, is a proinflammatory receptor. *J. Biol. Chem.* **288**(15), 10684-10691 (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.

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, 10/11/2022

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