

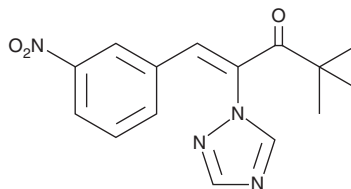
PRODUCT INFORMATION



Nexinhib20

Item No. 29899

CAS Registry No.: 331949-35-0
Formal Name: 4,4-dimethyl-1-(3-nitrophenyl)-2-(1H-1,2,4-triazol-1-yl)-1-penten-3-one
MF: $C_{15}H_{16}N_4O_3$
FW: 300.3
Purity: $\geq 98\%$
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

Nexinhib20 is supplied as a crystalline solid. A stock solution may be made by dissolving the nexinhib20 in the solvent of choice, which should be purged with an inert gas. Nexinhib20 is soluble in organic solvents such as ethanol and DMSO. The solubility of nexinhib20 in these solvents is approximately 15 and 30 mg/ml, respectively.

Description

Nexinhib20 is an inhibitor of the protein-protein interaction between the small GTPase Rab27a and its effector JFC1 ($IC_{50} = 2.6 \mu\text{M}$).¹ It inhibits myeloperoxidase (MPO) secretion induced by granulocyte macrophage colony-stimulating factor (GM-CSF) and fMLP in isolated human neutrophils ($IC_{50} = 0.33 \mu\text{M}$). Nexinhib20 (30 mg/kg) reduces LPS-induced increases in plasma MPO levels, as well as LPS-induced hepatic and renal neutrophil infiltration in a mouse model of endotoxin-induced systemic inflammation.

Reference

1. Johnson, J.L., Ramadass, M., He, J., *et al.* Identification of neutrophil exocytosis inhibitors (nexinhibs), small molecule inhibitors of neutrophil exocytosis and inflammation: Druggability of the small GTPase Rab27a. *J. Biol. Chem.* **291**(50), 25965-25982 (2016).

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, 12/09/2022

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