

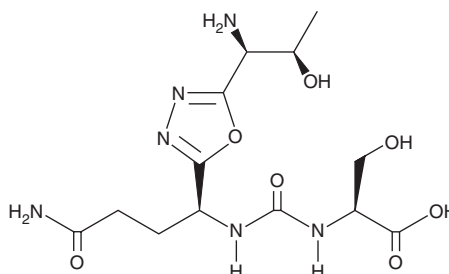
PRODUCT INFORMATION



PD-1-IN-17

Item No. 32797

CAS Registry No.: 1673560-66-1
Formal Name: N-[[[(1S)-4-amino-1-[5-[(1S,2R)-1-amino-2-hydroxypropyl]-1,3,4-oxadiazol-2-yl]-4-oxobutyl]amino]carbonyl]-L-serine
MF: C₁₃H₂₂N₆O₇
FW: 374.4
Purity: ≥95%
Supplied as: A 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

PD-1-IN-17 is supplied as a solid. A stock solution may be made by dissolving the PD-1-IN-17 in the solvent of choice, which should be purged with an inert gas. PD-1-IN-17 is soluble in the organic solvent DMSO.

Description

PD-1-IN-17 is an inhibitor of the programmed cell death protein 1 (PD-1) signaling pathway.¹ It reverses inhibition of isolated mouse splenocyte proliferation induced by PD-1 ligand (PD-L1) when used at a concentration of 100 nM.

Reference

1. Sasikumar, P.G.N., Ramachandra, M., and Naremaddepalli, S.S.S. 1,3,4-oxadiazole and 1,3,4-thiadiazole derivatives as immunomodulators. *Aurigen Discovery Technologies Limited*. **WO2015033301A1** (2015).

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, 10/17/2022

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