

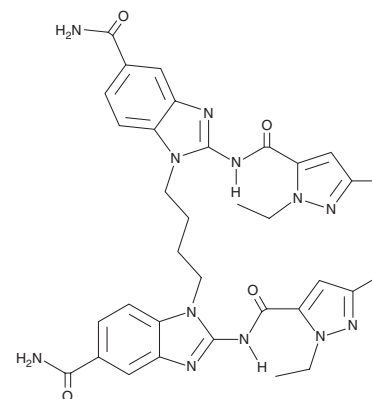
PRODUCT INFORMATION



diABZI STING Agonist 2

Item No. 34088

CAS Registry No.: 2138300-40-8
Formal Name: 1,1'-(1,4-butanediyl)bis[2-[[[(1-ethyl-3-methyl-1H-pyrazol-5-yl) carbonyl]amino]-1H-benzimidazole-5-carboxamide
Synonyms: CS-0087594, diABZI Stimulator of Interferon Genes Agonist 2, Diamidobenzimidazole STING Agonist-2, Diamidobenzimidazole Stimulator of Interferon Genes Agonist 2, HY-123943, STING Agonist diABZI Compound 2, STING Agonist (Compound 4), STING Agonist 4, Stimulator of Interferon Genes Agonist (Compound 4), Stimulator of Interferon Genes Agonist 4
MF: C₃₄H₃₈N₁₂O₄
FW: 678.7
Purity: ≥98%
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

diABZI STING agonist 2 is supplied as a solid. A stock solution may be made by dissolving the diABZI STING agonist 2 in the solvent of choice, which should be purged with an inert gas. diABZI STING agonist 2 is slightly soluble in DMSO.

Description

diABZI STING agonist 2 is an agonist of stimulator of interferon genes (STING).¹ It binds to STING with an apparent K_d value of 1.6 nM. It is selective for STING over the acute phase reactant orosomucoid 1 (ORM1; apparent K_d = 79 nM). diABZI STING agonist 2 increases phosphorylation of interferon regulatory factor 3 (IRF3) and STING in human peripheral blood mononuclear cells (PBMCs). It also induces secretion of IFN- β from PBMCs (EC_{50} = 3.1 μ M).

Reference

1. Ramanjulu, J.M., Pesiridis, G.S., Yang, J., *et al.* Design of amidobenzimidazole STING receptor agonists with systemic activity. *Nature* **564(7736)**, 439-443 (2018).

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, 07/19/2023

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