

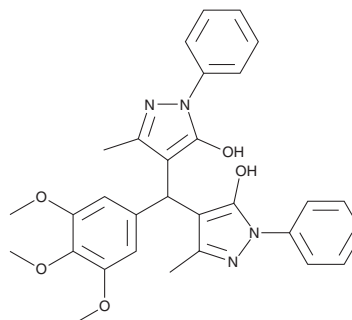
PRODUCT INFORMATION



SARS-CoV-2-IN-33

Item No. 45838

CAS Registry No.: 299919-79-2
Formal Name: 4,4'-[(3,4,5-trimethoxyphenyl)methylene]bis[3-methyl-1-phenyl-1H-pyrazol-5-ol]
Synonyms: ARS-CoV-2 Inhibitor 33, Severe Acute Respiratory Syndrome Coronavirus 2-IN-33, Severe Acute Respiratory Syndrome Coronavirus 2 Inhibitor 33
MF: C₃₀H₃₀N₄O₅
FW: 526.6
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

SARS-CoV-2-IN-33 is supplied as a solid. A stock solution may be made by dissolving the SARS-CoV-2-IN-33 in the solvent of choice, which should be purged with an inert gas. SARS-CoV-2-IN-33 is slightly soluble (0.1-1 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

Description

SARS-CoV-IN-33 is a pyrazolone derivative that has antiproliferative activity.¹ It inhibits the proliferation of MCF-7 and MDA-MB-231 breast, HeLa cervical, and PC3 prostate cancer cells (IC₅₀s = 5.45, 9.47, 22.86, and 28.44 μM, respectively) but not Ishikawa endometrial adenocarcinoma or HEK293 cells (IC₅₀s = >40 μM for both).

Reference

1. Gupta, A., Iqbal, S., Roohi, *et al.* Visible light-promoted green and sustainable approach for one pot synthesis of 4,4'-(arylmethylene)bis(1H-pyrazol-5-ols), in vitro anticancer activity, and molecular docking with Covid-19 M^{Pro}. *ACS Omega* **7**(38), 34583-34598 (2022).

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, 06/17/2026

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