

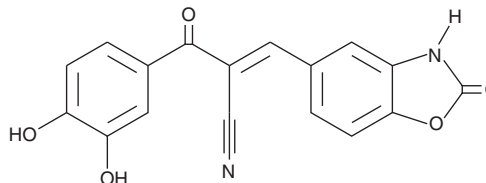
PRODUCT INFORMATION



AGL 2263

Item No. 21776

CAS Registry No.: 638213-98-6
Formal Name: α E-[(2,3-dihydro-2-oxo-5-benzoxazolyl)methylene]-3,4-dihydroxy- β -oxobenzenepropanenitrile
MF: C₁₇H₁₀N₂O₅
FW: 322.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 213, 222, 252, 350 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

AGL 2263 is supplied as a crystalline solid. A stock solution may be made by dissolving the AGL 2263 in the solvent of choice. AGL 2263 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of AGL 2263 in ethanol is approximately 2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

AGL 2263 is an inhibitor of the insulin-like growth factor 1 receptor (IGF-1R; IC₅₀ = 0.43 μ M).¹ It also inhibits the insulin receptor (IR), protein kinase B (PKB), and Src in a cell-free assay (IC₅₀s = 0.4, 55, and 2.2 μ M, respectively). AGL 2263 inhibits IGF-1R autophosphorylation and phosphorylation of the downstream elements IRS-1, PKB, and ERK2 in an ATP-independent manner. It inhibits colony formation of PC3 and LNCaP prostate and MCF-7 and MDA-MB-468 breast cancer cells in soft agar (IC₅₀s = 4.3, 9, 17, and 6 μ M, respectively).

Reference

1. Blum, G., Gazit, A., and Levitzki, A. Development of new insulin-like growth factor-1 receptor kinase inhibitors using catechol mimics. *J. Biol. Chem.* **278**(42), 40442-40454 (2003).

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, 12/22/2022

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