

PRODUCT INFORMATION



UNBS5162

Item No. 17739

CAS Registry No.: 956590-23-1

Formal Name: N-[2-[2-(dimethylamino)ethyl]-2,3-dihydro-1,3-dioxo-1H-benz[de]isoquinolin-5-yl]-urea

MF: C₁₇H₁₈N₄O₃

FW: 326.4

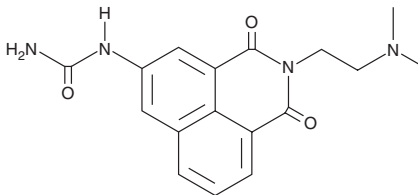
Purity: ≥98%

UV/Vis.: λ_{max}: 226, 249, 339, 384 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

UNBS5162 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNBS5162 in the solvent of choice, which should be purged with an inert gas. UNBS5162 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of UNBS5162 in these solvents is approximately 2 and 3 mg/ml, respectively.

UNBS5162 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UNBS5162 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. UNBS5162 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

UNBS5162 is a naphthalimide that antagonizes the expression of CXCL chemokines. *In vitro* exposure of PC-3 prostate cancer cells to 1 μM UNBS5162 for five successive days was shown to decrease the expression of proangiogenic CXCL chemokines.¹ UNBS5162 also demonstrates antiangiogenic properties *in vivo* in an orthotopic PC-3 model.¹

Reference

1. Mijatovic, T., Mahieu, T., Bruyère, C., *et al.* UNBS5162, a novel naphthalimide that decreases CXCL chemokine expression in experimental prostate cancers. *Neoplasia* **10**(6), 573-586 (2008).

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, 10/06/2022

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