

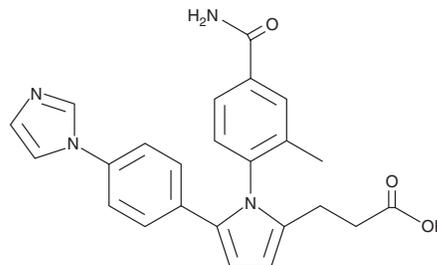
PRODUCT INFORMATION



N6022

Item No. 21269

CAS Registry No.: 1208315-24-5
Formal Name: 1-[4-(aminocarbonyl)-2-methylphenyl]-5-[4-(1H-imidazol-1-yl)phenyl]-1H-pyrrole-2-propanoic acid
MF: C₂₄H₂₂N₄O₃
FW: 414.5
Purity: ≥98%
UV/Vis.: λ_{max}: 304 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

N6022 is supplied as a crystalline solid. A stock solution may be made by dissolving the N6022 in the solvent of choice, which should be purged with an inert gas. N6022 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N6022 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of N6022 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of N6022 in PBS (pH 7.2) is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N6022 is a tight-binding, specific, and fully reversible inhibitor of S-nitrosogluthathione reductase (GSNOR; IC₅₀ = 8 nM; K_i = 2.5 nM).¹ It is reported to bind in the GSNO substrate binding pocket like a competitive inhibitor, although in kinetic assays it behaves with a mixed uncompetitive mode of inhibition toward the GSNO substrate and a mixed competitive mode of inhibition toward the formaldehyde adduct, S-hydroxymethylglutathione.¹

Reference

1. Green, L.S., Chun, L.E., Patton, A.K., *et al.* Mechanism of inhibition for N6022, a first-in-class drug targeting S-nitrosogluthathione reductase. *Biochemistry* **51(10)**, 2157-2168 (2012).

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, 12/07/2022

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