

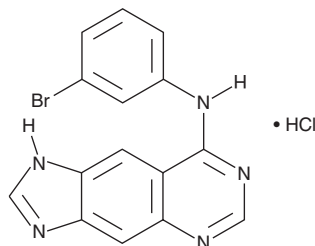
PRODUCT INFORMATION



BPIQ-II (hydrochloride)

Item No. 17226

CAS Registry No.: 171179-37-6
Formal Name: N-(3-bromophenyl)-3H-imidazo[4,5-g]quinazolin-8-amine, monohydrochloride
Synonym: PD 158294
MF: C₁₅H₁₀BrN₅ • HCl
FW: 376.6
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 349 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

BPIQ-II (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the BPIQ-II (hydrochloride) in the solvent of choice, which should be purged with an inert gas. BPIQ-II (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BPIQ-II (hydrochloride) in these solvents is approximately 30 and 10 mg/ml, respectively.

BPIQ-II (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BPIQ-II (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BPIQ-II (hydrochloride) has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BPIQ-II is a linear imidazoquinazoline that potently inhibits the tyrosine kinase activity of the epidermal growth factor receptor (EGFR; IC₅₀ = 8 pM).¹ It is selective for EGFR over an assortment of other tyrosine and serine/threonine kinases. Cellular studies indicate that BPIQ-II can enter cells and very selectively shut down EGF-stimulated signal transmission by binding competitively at the ATP site of EGFR.¹

Reference

1. Rewcastle, G.W., Palmer, B.D., Bridges, A.J., *et al.* Tyrosine kinase inhibitors. 9. Synthesis and evaluation of fused tricyclic quinazoline analogues as ATP site inhibitors of the tyrosine kinase activity of the epidermal growth factor receptor. *J. Med. Chem.* **39**(4), 918-928 (1996).

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, 12/14/2022

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