

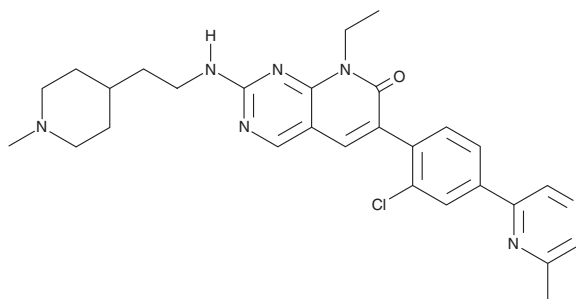
PRODUCT INFORMATION



FRAX1036

Item No. 20711

CAS Registry No.: 1432908-05-8
Formal Name: 6-[2-chloro-4-(6-methyl-2-pyrazinyl)phenyl]-8-ethyl-2-[[2-(1-methyl-4-piperidyl)ethyl]amino]-pyrido[2,3-d]pyrimidin-7(8H)-one
MF: C₂₈H₃₂ClN₇O
FW: 518.1
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 299, 308, 355 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

FRAX1036 is supplied as a crystalline solid. A stock solution may be made by dissolving the FRAX1036 in the solvent of choice, which should be purged with an inert gas. FRAX1036 is soluble in organic solvents such as ethanol and dimethyl formamide (DMF). The solubility of FRAX1036 in these solvents is approximately 0.33 and 1 mg/ml, respectively.

FRAX1036 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, FRAX1036 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. FRAX1036 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

FRAX1036 is a selective inhibitor of p21-activated kinase 1 (PAK1), a serine/threonine kinase downstream of Rac1 and Cdc42 that is involved in tumorigenesis. It can induce apoptosis in breast cancer cells and has been shown to potentiate the activity of the microtubule stabilizing agent, docetaxel (Item No. 11637).¹ Disruption of PAK1 via FRAX1036 has been used to inhibit oncogenic KRAS signaling in non-small cell lung cancer cells.²

References

1. Ong, C.C., Gierke, S., Pitt, C., *et al.* Small molecule inhibition of group I p21-activated kinases in breast cancer induces apoptosis and potentiates the activity of microtubule stabilizing agents. *Breast Cancer Res.* **17(59)** (2015).
2. Mortazavi, F., Lu, J., Phan, R., *et al.* Significance of KRAS/PAK1/Crk pathway in non-small cell lung cancer oncogenesis. *BMC Cancer* **15(381)** (2015).

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, 12/21/2022

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