

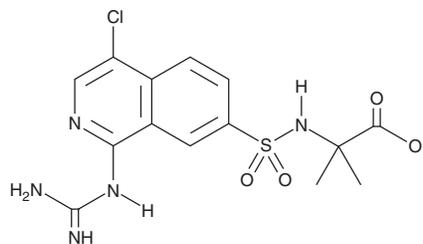
PRODUCT INFORMATION



UK 371804

Item No. 27225

CAS Registry No.: 256477-09-5
Formal Name: N-[[1-[(aminoiminomethyl)amino]-4-chloro-7-isoquinolinyl]sulfonyl]-2-methyl-alanine
MF: C₁₄H₁₆ClN₅O₄S
FW: 385.8
Purity: ≥95%
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

UK 371804 is supplied as a crystalline solid. A stock solution may be made by dissolving the UK 371804 in the solvent of choice, which should be purged with an inert gas. UK 371804 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

UK 371804 is an inhibitor of urokinase-type plasminogen activator (uPA; $K_i = 10 \text{ nM}$).¹ It is 4,000- and 2,700-fold selective for uPA over tissue plasminogen activator (tPA) and plasmin, respectively. UK 371804 inhibits exogenous uPA in human chronic wound fluid ($IC_{50} = 0.89 \text{ } \mu\text{M}$). Topical administration of UK 371804 inhibits exogenous uPA activity in a porcine excisional wound model.

Reference

1. Fish, P.V., Barber, C.G., Brown, D.G., *et al.* Selective urokinase-type plasminogen activator inhibitors. 4. 1-(7-Sulfonamidoisoquinolinyl)guanidines. *J. Med. Chem.* **50**(10), 2341-2351 (2007).

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, 11/09/2022

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