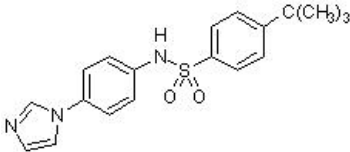


Product datasheet

ISCK03, SCF/c-Kit inhibitor ab145066

1 Image

Overview

Product name	ISCK03, SCF/c-Kit inhibitor
Description	Cell-permeable, selective SCF/c-Kit inhibitor
Biological description	Cell-permeable, selective SCF/c-Kit inhibitor (IC ₅₀ < 2.5 μM). Blocks SCF-induced c-kit and Erk phosphorylation without affecting HGF-induced Erk phosphorylation. Shows antiangiogenic effects. Abolishes melanin production <i>in vivo</i> .
Purity	> 99%
CAS Number	945526-43-2
Chemical structure	

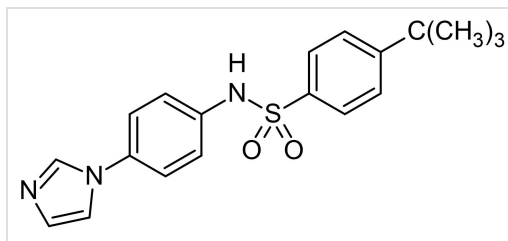
Properties

Chemical name	4-(1,1-Dimethylethyl-N-[4-(1H-imidazol-1-yl)phenyl]benzenesulfonamide
Molecular weight	355.45
Molecular formula	C ₁₉ H ₂₁ N ₃ O ₂ S
PubChem identifier	3792751
Storage instructions	Store at +4°C. Store under desiccating conditions. The product can be stored for up to 12 months.
Solubility overview	Soluble in ethanol to 25 mM and in DMSO to 100 mM
Handling	Wherever possible, you should prepare and use solutions on the same day. However, if you need to make up stock solutions in advance, we recommend that you store the solution as aliquots in tightly sealed vials at -20°C. Generally, these will be useable for up to one month. Before use, and prior to opening the vial we recommend that you allow your product to equilibrate to room temperature for at least 1 hour. Need more advice on solubility, usage and handling? Please visit our frequently asked questions (FAQ) page for more details.
SMILES	CC(C)(C)C1=CC=C(C=C1)S(=O)(=O)NC2=CC=C(C=C2)N3C=CN=C3

Source

Synthetic

Images



2D chemical structure image of ab145066, ISCK03, SCF/c-Kit inhibitor

Chemical Structure - ISCK03, SCF/c-Kit inhibitor
(ab145066)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES, NOT FOR USE IN HUMANS"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors
- Abcam biochemicals are novel compounds and we have not tested their biological activity in house. Please use the literature to identify how to use these products effectively. If you require further assistance please contact the scientific support team