

Regorafenib

Multi-kinase inhibitor (Including KDR-Tie2)

Catalog # K01-900

Lot # D3905-3

CAS # 755037-03-7

Product Description

Molecular Formula: C₂₁H₁₅ClF₄N₄O₃

Appearance: Off-White powder

Molecular Weight: 482.80

Purity: >98% (HPLC,TLC); NMR (conforms)

Solubilization: May be dissolved in DMSO (50 mg/ml)

Alias

4-[4-[[[4-chloro-3-(trifluoromethyl)phenyl]amino]carbonyl]amino]-3-fluorophenoxy]-N-methyl-2-pyridinecarboxamide; BAY 73-4506

Specific Activity

Multi-kinase inhibitor (Including KDR-Tie2)

Storage and Stability

Store desiccated as supplied at -20°C for up to 2 year.

Store solutions at -20°C for up to 3 months.

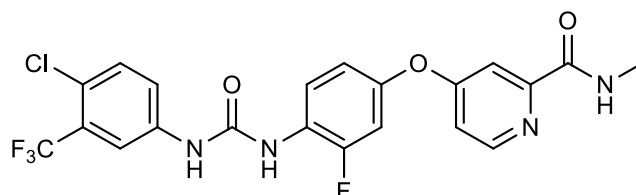
Scientific Background

A multi-kinase inhibitor related to sorafenib¹. Inhibits c-Raf, B-Raf, VEGFR1, VEGFR2, VEGFR3, Tie2, PDGFRβ, FGFR1, c-Kit, and RET with IC₅₀s=2.5, 28, 13, 4.2, 46, 311, 22, 202, 7 and 1.5 nM respectively¹⁻²

References

1. SM Wilhelm *et al.* *Int. J. Cancer* 2011 129:245
2. JCM Uitdehaag *et al.* *PLoS One* 2014 9(3):e92146

Molecular Structure



Regorafenib

Multi-kinase inhibitor (Including KDR-Tie2)

Catalog #	K01-900
CAS #	755037-03-7
Lot #	D3905-3
Purity	>98% (HPLC,TLC); NMR (conforms)
Format	Off-White powder
Solubilization	May be dissolved in DMSO (50 mg/ml)
Storage & Shipping	Store desiccated as supplied at -20°C for up to 1 years. Store solutions at -20°C for up to 3 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.