

# BAY 43-9006

Broad-spectrum kinase inhibitor

Broad-spectrum kinase inhibitor.

Citations: 5

[View Online »](#)

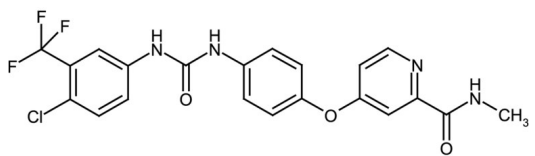
## Ordering Information

[Order Online »](#)

ALX-270-469-M001	1mg
ALX-270-469-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** N-[4-Chloro-3-(trifluoromethyl)phenyl]-([4-[2-(N-methylcarbamoyl)(4-pyridyloxy)]phenyl]amino)-carboxamide, Sorafenib

**Appearance** White to off-white solid.

**CAS** 284461-73-0

**Couple Target** Kinase

**Couple Type** Inhibitor

**Formula**  $C_{21}H_{16}ClF_3N_4O_3$

**Identity** Identity determined by 1H-NMR.

**MW** 464.8

**Purity** ≥95% (HPLC)

**Solubility** Soluble in DMSO, 100% ethanol, methanol or ethyl acetate.

**Source** Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)