

# PRODUCT INFORMATION



## AC-710

Item No. 29498

**CAS Registry No.:** 1351522-04-7  
**Formal Name:** N-[4-[[[5-(1,1-dimethylethyl)-3-isoxazolyl]amino]carbonyl]amino]phenyl]-5-[(1-ethyl-2,2,6,6-tetramethyl-4-piperidinyloxy)-2-pyridinecarboxamide

**MF:** C<sub>31</sub>H<sub>42</sub>N<sub>6</sub>O<sub>4</sub>  
**FW:** 562.7

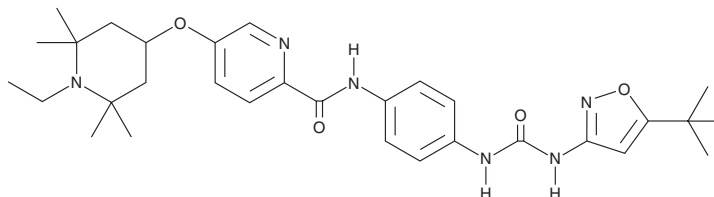
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 255, 299 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

AC-710 is supplied as a crystalline solid. A stock solution may be made by dissolving the AC-710 in the solvent of choice, which should be purged with an inert gas. AC-710 is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

### Description

AC-710 is an inhibitor of PDGFR family kinases (IC<sub>50</sub>s = 7.7, 10.5, 2, and 1.2 nM for PDGFRβ, CSF1R, FLT3, and c-Kit, respectively).<sup>1</sup> It is greater than 30-fold selective for PDGFR family kinases over a panel of 386 kinases, as well as over a panel of five cytochrome P450 (CYP) enzymes (IC<sub>50</sub>s = >40 μM). AC-710 (30 mg/kg) inhibits CSF1R-dependent M-NFS-60 murine leukemia cell proliferation and reduces tumor volume in an FLT3 mutant MV4-11 leukemia mouse xenograft model.

### Reference

1. Liu, G., Campbell, B.T., Holladay, M.W., *et al.* Discovery of AC710, a globally selective inhibitor of platelet-derived growth factor receptor-family kinases. *ACS Med. Chem. Lett.* **3**(12), 997-1002 (2012).

#### 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, 10/25/2022

#### 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