

# PRODUCT INFORMATION

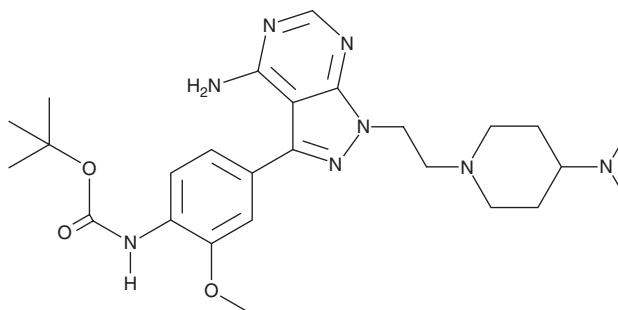


## eCF506

Item No. 19959

**CAS Registry No.:** 1914078-41-3  
**Formal Name:** N-[4-[4-amino-1-[2-[4-(dimethylamino)-1-piperidinyl]ethyl]-1H-pyrazolo[3,4-d]pyrimidin-3-yl]-2-methoxyphenyl]-carbamic acid, 1,1-dimethylethyl ester

**MF:** C<sub>26</sub>H<sub>38</sub>N<sub>8</sub>O<sub>3</sub>  
**FW:** 510.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 251, 268, 300 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

eCF506 is supplied as a crystalline solid. A stock solution may be made by dissolving the eCF506 in the solvent of choice, which should be purged with an inert gas. eCF506 is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of eCF506 in these solvents is approximately 15 mg/ml.

### Description

eCF506 is an inhibitor of Src kinases (IC<sub>50</sub>s = <0.5, 2.1, <0.5 nM for Src, Fyn, and Yes, respectively).<sup>1</sup> It is selective for Src kinases over Abl (IC<sub>50</sub> = 479 nM), as well as KIT, mTOR, PDGFRα, and RET (IC<sub>50</sub>s = >100 μM for all). eCF506 completely inhibits phosphorylation of Src and FAK, a downstream Src substrate, in MCF-7 and MDA-MB-231 cells when used at a concentration of 100 nM. It also inhibits proliferation of MCF-7 cells (EC<sub>50</sub> = 9 nM) and induces apoptosis. eCF506 (50 mg/kg per day) reduces levels of activated Src kinase in tumor tissue in an HCT116 mouse xenograft model.

### Reference

1. Fraser, C., Dawson, J.C., Dowling, R., *et al.* Rapid discovery and structure-activity relationships of pyrazolopyrimidines that potently suppress breast cancer cell growth via SRC kinase inhibition with exceptional selectivity over ABL kinase. *J. Med. Chem.* **59**(10), 4697-4710 (2016).

#### 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, 12/19/2022

#### 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](http://WWW.CAYMANCHEM.COM)