

# PRODUCT INFORMATION



## Bayer 18

Item No. 41124

CAS Registry No.: 1251752-12-1

Formal Name: N-(1,1-dimethylethyl)-N'-[3-[[5-fluoro-4-[(2-hydroxy-1,1-dimethylethyl)amino]-2-pyrimidinyl]amino]phenyl]-urea

MF:  $C_{19}H_{27}FN_6O_2$

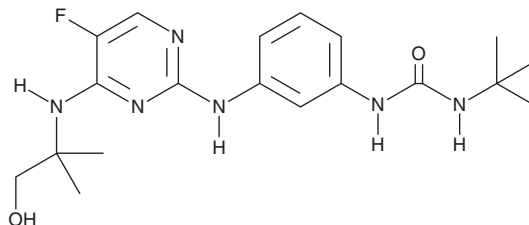
FW: 390.5

Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Bayer 18 is supplied as a solid. A stock solution may be made by dissolving the bayer 18 in the solvent of choice, which should be purged with an inert gas. Bayer 18 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

### Description

Bayer 18 is a tyrosine kinase 2 (TYK2) inhibitor ( $\text{IC}_{50} = 18.7 \text{ nM}$ ).<sup>1</sup> It is selective for TYK2 over JAK2, cyclin-dependent kinase 2 (Cdk2), and VEGFR2 ( $\text{IC}_{50}\text{s} = 2,000, 1,790, \text{ and } 3,520 \text{ nM}$ , respectively).

### Reference

1. Buchmann, B., Eis, K., Bothe, U., *et al.* Aminocabonylamino-substituted anilino-pyrimidine derivatives as Tyk inhibitors, their production and use as medicinal products. **DE 10 2009 015 070 A1** (2010).

#### 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, 07/09/2024

#### 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