

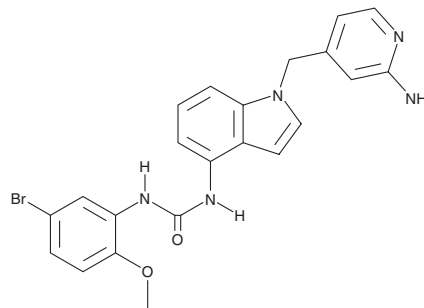
# PRODUCT INFORMATION



## Jl-101

Item No. 36405

**CAS Registry No.:** 900573-88-8  
**Formal Name:** N-[1-[(2-amino-4-pyridinyl)methyl]-1H-indol-4-yl]-N'-(5-bromo-2-methoxyphenyl)-urea  
**Synonym:** CGI1842  
**MF:** C<sub>22</sub>H<sub>20</sub>BrN<sub>5</sub>O<sub>2</sub>  
**FW:** 466.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 221, 308 nm  
**Supplied as:** A 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

Jl-101 is supplied as a solid. A stock solution may be made by dissolving the Jl-101 in the solvent of choice, which should be purged with an inert gas. Jl-101 is soluble in DMSO and methanol (heated).

### Description

Jl-101 is an inhibitor of VEGFR2, PDGFRβ, and Eph tyrosine kinase receptor 4 (EphB4) with anticancer activity.<sup>1</sup>

### Reference

1. Velleca, M., Miller, L., Sharma, S., *et al.* Abstract 17: Preclinical and phase 1 trial results of Jl-101, a novel, oral tyrosine kinase inhibitor that selectively targets VEGFR2, EphB4, and PDGFRβ. *Cancer Res.* 70(Suppl. 8), 17 (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, 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