

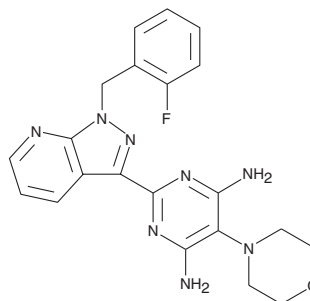
# PRODUCT INFORMATION



## BAY 41-8543

Item No. 10011131

**CAS Registry No.:** 256498-66-5  
**Formal Name:** 2-[1-[(2-fluorophenyl)methyl]-1H-pyrazolo[3,4-b]pyridin-3-yl]-5-(4-morpholinyl)-4,6-pyrimidinediamine  
**MF:** C<sub>21</sub>H<sub>21</sub>FN<sub>8</sub>O  
**FW:** 420.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212, 327 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

BAY 41-8543 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY 41-8543 in the solvent of choice, which should be purged with an inert gas. BAY 41-8543 is soluble in dimethyl formamide (DMF). The solubility of BAY 41-8543 in this solvent is approximately 3 mg/ml.

BAY 41-8543 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAY 41-8543 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BAY 41-8543 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Soluble guanylate cyclase (sGC) is the primary cellular receptor for nitric oxide (NO). NO binds and activates a heme group in sGC, initiating the conversion of GTP to the second messenger cyclic GMP (cGMP). BAY 41-8543 is a heme-dependent stimulator of sGC, increasing the activity of recombinant sGC dose-dependently, from 0.1 nM to 100 μM, up to 92-fold.<sup>1</sup> Surprisingly, NO donors synergize with BAY 41-8543 in stimulating recombinant sGC.<sup>1</sup> BAY 41-8543 relaxes vessels and inhibits platelet aggregation *in vitro* at nM concentrations.<sup>1</sup> *In vivo*, BAY 41-8543 decreases blood pressure dose-dependently, prolongs bleeding time, and reduces thrombosis.<sup>2</sup> Inhalation of microparticles containing BAY 41-8543 increases pulmonary vasodilation without changing mean arterial pressure, suggesting that agonists of sGC may be efficacious in treating pulmonary hypertension.<sup>3</sup>

### References

1. Stasch, J.-P., Alonso-Alija, C., Apeler, H., *et al.* Pharmacological actions of a novel NO-independent guanylyl cyclase stimulator, BAY 41-8543: *In vitro* studies. *Br. J. Pharmacol.* **135**(2), 333-343 (2002).
2. Stasch, J.-P., Dembowski, K., Perzborn, E., *et al.* Cardiovascular actions of a novel NO-independent guanylyl cyclase stimulator, BAY 41-8543: *In vivo* studies. *Br. J. Pharmacol.* **135**(2), 344-355 (2002).
3. Evgenov, O.V., Kohane, D.S., Bloch, K.D., *et al.* Inhaled agonists of soluble guanylate cyclase induce selective pulmonary vasodilation. *Am. J. Respir. Crit. Care Med.* **176**(11), 1138-1145 (2007).

**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, 01/31/2024

### 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