

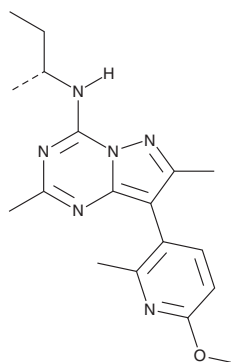
# PRODUCT INFORMATION



## Pexacerfont

Item No. 33221

**CAS Registry No.:** 459856-18-9  
**Formal Name:** 8-(6-methoxy-2-methyl-3-pyridinyl)-2,7-dimethyl-N-[(1R)-1-methylpropyl]-pyrazolo[1,5-a]-1,3,5-triazin-4-amine  
**Synonyms:** BMS 562086, DPC-A69448  
**MF:** C<sub>18</sub>H<sub>24</sub>N<sub>6</sub>O  
**FW:** 340.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 283 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

Pexacerfont is supplied as a crystalline solid. A stock solution may be made by dissolving the pexacerfont in the solvent of choice, which should be purged with an inert gas. Pexacerfont is soluble in organic solvents such as ethanol and DMSO. The solubility of pexacerfont in these solvents is approximately 30 mg/ml.

### Description

Pexacerfont is an antagonist of corticotropin-releasing factor receptor 1 (CRF<sub>1</sub>; IC<sub>50</sub> = 6.1 nM for the human receptor).<sup>1</sup> It is selective for CRF<sub>1</sub> in IMR-32 cells (mean IC<sub>50</sub> = 10.6 nM) over CFR<sub>2β</sub> receptors endogenously expressed in pig choroid plexus membranes (IC<sub>50</sub> = >1,000 nM) in radioligand binding assays and does not bind to the human CRF binding protein (IC<sub>50</sub> = >1,000 nM). Pexacerfont inhibits adrenocorticotrophic hormone (ACTH) release induced by CRF in primary rat pituitary cells with an IC<sub>50</sub> value of 129 nM. It increases the time spent in the open arms of the elevated plus maze (EPM) in rats, indicating anxiolytic-like activity, when administered at doses ranging from 10 to 30 mg/kg.

### Reference

1. Gilligan, P.J., Clarke, T., He, L., *et al.* Synthesis and structure-activity relationships of 8-(pyrid-3-yl)pyrazolo[1,5-a]-1,3,5-triazines: Potent, orally bioavailable corticotropin releasing factor receptor-1 (CRF<sub>1</sub>) antagonists. *J. Med. Chem.* **52(9)**, 3084-3092 (2009).

#### 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, 11/14/2022

#### 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