

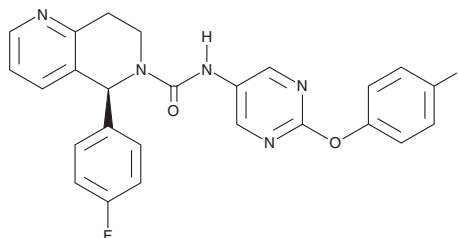
# PRODUCT INFORMATION



## BAY-899

Item No. 41716

**CAS Registry No.:** 2471967-92-5  
**Formal Name:** N-[2-(4-fluorophenoxy)-5-pyrimidinyl]-5-(4-fluorophenyl)-7,8-dihydro-1,6-naphthyridine-6(5H)-carboxamide  
**MF:** C<sub>25</sub>H<sub>19</sub>F<sub>2</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 459.5  
**Purity:** ≥98%  
**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

BAY-899 is supplied as a solid. A stock solution may be made by dissolving the BAY-899 in the solvent of choice, which should be purged with an inert gas. BAY-899 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

BAY-899 is an antagonist of the luteinizing hormone receptor (LHR; IC<sub>50</sub>s = 185 and 46 nM for the human and rat receptors, respectively).<sup>1</sup> It is selective for LHR over thyroid-stimulating hormone (TSH) and follicle-stimulating hormone (FSH) receptors (IC<sub>50</sub>s = 24 and >16 μM, respectively, for the human receptors) and a panel of 25 G protein-coupled receptors (GPCRs) at 10 μM. BAY-899 (12.5 mg/kg per day) reduces serum levels of estradiol in female rats.

### Reference

1. Wortmann, L., Lindenthal, B., Muhn, P., *et al.* Discovery of BAY-298 and BAY-899: Tetrahydro-1,6-naphthyridine-based, potent, and selective antagonists of the luteinizing hormone receptor which reduce sex hormone levels in vivo. *J. Med. Chem.* **62**(22), 10321-10341 (2019).

#### 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, 09/06/2024

#### 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