

PRODUCT INFORMATION



YIL 781 (hydrochloride)

Item No. 29217

CAS Registry No.: 1640226-17-0

Formal Name: 6-(4-fluorophenoxy)-2-methyl-3-[[[(3S)-1-(1-methylethyl)-3-piperidiny]methyl]-4(3H)-quinazolinone, monohydrochloride

MF: $C_{24}H_{28}FN_3O_2 \cdot HCl$

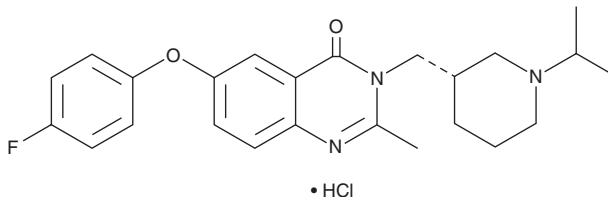
FW: 446.0

Purity: $\geq 98\%$

Supplied as: A solid

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

Stability: ≥ 4 years



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

Laboratory Procedures

YIL 781 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the YIL 781 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. YIL 781 (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. The solubility of YIL 781 (hydrochloride) in these solvents is approximately 100 mM.

Description

YIL 781 is an antagonist of the growth hormone secretagogue receptor 1a (GHS-R1a; $K_i = 0.017 \mu M$).¹ It is selective for GHS-R1a over the motilin receptor ($K_i = 6 \mu M$) and a panel of 30 receptors and enzymes (K_i s = $>10 \mu M$). YIL 781 (1 μM) prevents ghrelin-induced decreases in insulin secretion in glucose-stimulated isolated rat pancreatic islets. It increases plasma insulin levels and decreases blood glucose levels in an intraperitoneal glucose tolerance test in rats when administered at doses of 0.3 or 10 mg/kg, respectively. YIL 781 (10 and 20 mg/kg) inhibits cocaine-induced hyperlocomotion in mice.²

References

1. Esler, W.P., Rudolph, J., Claus, T.H., *et al.* Small-molecule ghrelin receptor antagonists improve glucose tolerance, suppress appetite, and promote weight loss. *Endocrinology* **148**(11), 5175-5185 (2007).
2. Toth, K., Slosky, L.M., Pack, T.F., *et al.* Ghrelin receptor antagonism of hyperlocomotion in cocaine-sensitized mice requires β arrestin-2. *Synapse* **72**(1), e22012 (2018).

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, 10/28/2022

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