

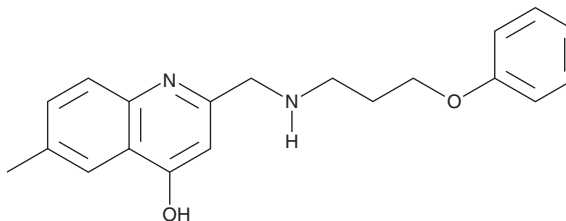
# PRODUCT INFORMATION



## UCSF924

Item No. 41317

CAS Registry No.: 1434515-70-4  
Formal Name: 6-methyl-2-[[[3-phenoxypropyl]amino]methyl]-4-quinolinol  
MF:  $C_{20}H_{22}N_2O_2$   
FW: 322.4  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

UCSF924 is supplied as a solid. A stock solution may be made by dissolving the UCSF924 in the solvent of choice, which should be purged with an inert gas. UCSF924 is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

### Description

UCSF924 is a dopamine  $D_4$  receptor agonist.<sup>1</sup> It selectively induces  $\beta$ -arrestin translocation over  $G_{\alpha_{i/o}}$  signaling in HEK293T cells expressing human dopamine  $D_4$  receptors by 7.4-fold at 1  $\mu\text{M}$ . UCSF924 selectively induces cAMP accumulation in HEK293T cells expressing human dopamine  $D_4$  receptors ( $EC_{50} = 4.2$  nM) over HEK293T cells expressing human dopamine  $D_2$  or dopamine  $D_3$  receptors ( $EC_{50}$ s =  $>10$   $\mu\text{M}$  for both) and induces  $\beta$ -arrestin recruitment in HTLA cells expressing human dopamine  $D_4$  receptors over a panel of HTLA cells expressing one of 320 non-olfactory G protein-coupled receptors (GPCRs) at 1  $\mu\text{M}$ .

### Reference

1. Wang, S., Wacker, D., Levit, A., et al.  $D_4$  dopamine receptor high-resolution structures enable the discovery of selective agonists. *Science* **358**(6361), 381-386 (2017).

#### 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, 07/01/2024

#### 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