

# PRODUCT INFORMATION



## L-742,001 (hydrochloride)

Item No. 40920

CAS Registry No.: 174605-64-2

Formal Name: 4-[4-[(4-chlorophenyl)methyl]-1-(phenylmethyl)-4-piperidinyl]-2-hydroxy-4-oxo-2-butenoic acid, monohydrochloride

MF:  $C_{23}H_{24}ClNO_4 \cdot HCl$

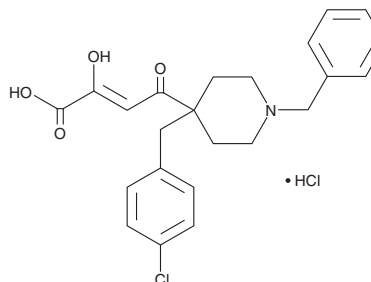
FW: 450.4

Purity:  $\geq 98\%$

Supplied as: A solid

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

Stability:  $\geq 4$  years



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

### Laboratory Procedures

L-742,001 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the L-742,001 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. L-742,001 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

### Description

L-742,001 is an inhibitor of influenza virus PA endonuclease ( $IC_{50} = 0.5 \mu M$ ), an essential subdomain of the viral RNA polymerase.<sup>1</sup> It reduces viral titers in PR8 influenza-infected MDCK cells when used at a concentration of  $10 \mu M$ .<sup>2</sup>

### References

1. Stevaert, A., Nurra, S., Pala, N., *et al.* An integrated biological approach to guide the development of metal-chelating inhibitors of influenza virus PA endonuclease. *Mol. Pharmacol.* **87**(2), 323-337 (2015).
2. Song, M.-S., Kumar, G., Shadrack, W.R., *et al.* Identification and characterization of influenza variants resistant to a viral endonuclease inhibitor. *Proc. Nat. Acad. Sci. USA* **113**(13), 3669-3674 (2016).

#### 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, 05/17/2024

#### 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