PRODUCT INFORMATION



Gisadenafil (besylate)

Item No. 31472

CAS Registry No.:	334827-98-4	0
Formal Name:	5-[2-ethoxy-5-[(4-ethyl-	
	1-piperazinyl)sulfonyl]-3-	
	pyridinyl]-3-ethyl-2,6-dihydro-	
	2-(2-methoxyethyl)-7H-	
	pyrazolo[4,3-d]pyrimidin-7-one,	
	monobenzenesulfonate	
Synonym:	UK369003	
MF:	$C_{23}H_{33}N_7O_5S \bullet C_6H_6O_3S$	0_0
FW:	677.8	Ś
Purity:	≥98%	. [\ OH
Supplied as:	A solid	
Storage:	-20°C	\checkmark
Stability:	≥4 years	

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

Laboratory Procedures

Gisadenafil (besylate) is supplied as a solid. A stock solution may be made by dissolving the gisadenafil (besylate) in the solvent of choice, which should be purged with an inert gas. Gisadenafil (besylate) is soluble in the organic solvent DMSO. It is also soluble in water. The solubility of gisadenafil (besylate) in DMSO and water is approximately 50 and 5 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Gisadenafil is a phosphodiesterase 5 (PDE5) inhibitor (IC₅₀ = 1.1 nM).¹ It is 103-fold selective for PDE5 over PDE6 at 1.1 nM. Gisadenafil (2 mg/kg) restores the wild-type response to hypercapnia, increased cortical blood flow and dilation of small arterioles, in the GFAP-driven, doxycycline-inducible HIV-1 Tat transgenic (Tat-tg) mouse model of HIV-1 neuroinflammation.²

References

- 1. Rawson, D.J., Ballard, S., Barber, C., et al. The discovery of UK-369003, a novel PDE5 inhibitor with the potential for oral bioavailability and dose-proportional pharmacokinetics. Bioorg. Med. Chem. 20(1), 498-509 (2012).
- 2. Silva, J., Polesskaya, O., Knight, W., et al. Transient hypercapnia reveals an underlying cerebrovascular pathology in a murine model for HIV-1 associated neuroinflammation: role of NO-cGMP signaling and normalization by inhibition of cyclic nucleotide phosphodiesterase-5. J. Neuroinflammation 9, 253 (2012).

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.

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, 11/10/2022

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