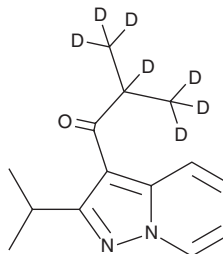


PRODUCT INFORMATION



Ibudilast-d₇ Item No. 28876

CAS Registry No.: 2713301-45-0
Formal Name: 2-(methyl-d₃)-1-[2-(1-methylethyl)pyrazolo[1,5-a]pyridin-3-yl]-1-propanone-2,3,3,3-d₄
Synonyms: AV 411-d₇, KC-404-d₇
MF: C₁₄H₁₁D₇N₂O
FW: 237.4
Chemical Purity: ≥98% (Ibudilast)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
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

Ibudilast-d₇ is intended for use as an internal standard for the quantification of ibudilast (Item No. 14832) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Ibudilast-d₇ is supplied as a solid. A stock solution may be made by dissolving the ibudilast-d₇ in the solvent of choice, which should be purged with an inert gas. Ibudilast-d₇ is soluble in chloroform and methanol.

Description

Ibudilast is an inhibitor of phosphodiesterase 4 (PDE4; IC₅₀s = 54, 65, 239, and 166 nM for PDE4A-D, respectively).¹ It is selective for PDE4 over PDE1, PDE7A, PDE7B, and PDE9A (IC₅₀s = ≥10,000 nM for all) but does inhibit PDE3A, PDE3B, and PDE5A (IC₅₀s = 1,600, 2,700, and 3,510 nM, respectively). Ibudilast inhibits LPS-induced production of TNF-α and fMLP-induced production of leukotriene B₄ (LTB₄; Item No. 20110) in isolated human whole blood (IC₅₀s = 6.2 and 2.5 μM, respectively). It inhibits bronchospasm by 34% in a guinea pig model of leukotriene-mediated allergic bronchospasm when administered intravenously at a dose of 5 mg/kg.² Ibudilast prevents increases in TNF-α, IL-1β, and IL-6 expression in the striatum in a mouse model of MPTP-induced Parkinson's disease.³ It also increases striatal expression of glial cell-derived neurotrophic factor (GDNF) in MPTP-treated and -untreated mice when administered at doses of 40 and 50 mg/kg, respectively, twice per day.

References

1. Huang, Z., Liu, S., Zhang, L., *et al.* Preferential inhibition of human phosphodiesterase 4 by ibudilast. *Life Sci.* **78(23)**, 2663-2668 (2006).
2. Kreutner, W., Sherwood, J., and Rizzo, C. The effect of leukotriene antagonists, lipoxygenase inhibitors and selected standards on leukotriene-mediated allergic bronchospasm in guinea pigs. *Agents Actions* **28(3-4)**, 173-184 (1989).
3. Schwenkgrub, J., Zaremba, M., Joniec-Maciejak, I., *et al.* The phosphodiesterase inhibitor, ibudilast, attenuates neuroinflammation in the MPTP model of Parkinson's disease. *PLoS One* **12(7)**, e0182019 (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, 03/08/2023

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