

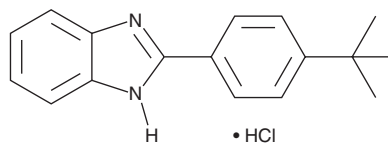
PRODUCT INFORMATION



ZLN005 (hydrochloride)

Item No. 9002449

Formal Name: 2-[4-(1,1-dimethylethyl)phenyl]-1H-benzimidazole, monohydrochloride
MF: C₁₇H₁₈N₂ • HCl
FW: 286.8
Purity: ≥98%
UV/Vis.: λ_{max}: 250, 303 nm
Supplied as: A crystalline 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

ZLN005 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the ZLN005 (hydrochloride) in the solvent of choice. ZLN005 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of ZLN005 (hydrochloride) in these solvents is approximately 25, 17, and 11 mg/ml, respectively.

ZLN005 (hydrochloride) is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Peroxisome proliferator-activated receptor-γ coactivator-1α (PGC-1α) is a tissue-specific and inducible transcriptional coactivator for several nuclear receptors and plays a key role in energy metabolism, hepatic gluconeogenesis, and cholesterol homeostasis.¹ ZLN005 is a small molecule that stimulates the expression of PGC-1α and downstream genes in skeletal muscle cells, improving glucose utilization and fatty acid oxidation at a concentration of 20 μM.² Chronic administration of 15 mg/kg/day ZLN005 to diabetic *db/db* mice increased PGC-1α and downstream gene transcription in skeletal muscle, increasing fat oxidation and improving glucose tolerance, pyruvate tolerance, and insulin sensitivity.² This compound is the hydrochloride salt formulation of Item ZLN005 (Item No. 14121).

References

1. Lin, J., Handschin, C., and Spiegelman, B.M. Metabolic control through the PGC-1 family of transcription coactivators. *Cell Metab.* **1**, 361-370 (2005).
2. Zhang, L.-N., Zhou, H.-Y., Fu, Y.-Y., et al. Novel small-molecule PGC-1α transcriptional regulator with beneficial effects on diabetic *db/db* mice. *Diabetes* **62**(4), 1297-1307 (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.

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, 05/01/2024

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