

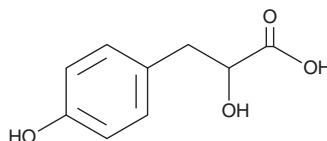
PRODUCT INFORMATION



3-(4-Hydroxyphenyl)lactic Acid

Item No. 36367

CAS Registry No.: 306-23-0
Formal Name: α ,4-dihydroxy-benzenepropanoic acid
Synonyms: HPLA, ($\pm$)-3-(4-Hydroxyphenyl)lactic Acid,
 β -(4-Hydroxyphenyl)-DL-Lactic Acid,
DL-*p*-Hydroxyphenyllactic Acid
MF: $C_9H_{10}O_4$
FW: 182.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 226 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

3-(4-Hydroxyphenyl)lactic acid (HPLA) is supplied as a solid. A stock solution may be made by dissolving the HPLA in the solvent of choice, which should be purged with an inert gas. HPLA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of HPLA in these solvents is approximately 25 mg/ml.

Description

HPLA is a bacterial metabolite produced by *L. murinus* in the gut microbiome.¹ Oral administration of HPLA with DL-3-phenylacetic acid (PLA), 2-hydroxyisocaproic acid (HICA), 2-hydroxy-3-methylbutyric acid (HMBA), and 1-indole-lactic acid (ILA) increases intestinal and white adipose tissue lipid absorption and decreases fatty acid oxidation in a mouse model of refeeding-induced obesity.

Reference

1. Zarrinpar, A. A high-protein diet prevents weight regain. *Nat. Metab.* **4**(12), 1616-1617 (2022).

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, 02/06/2023

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