

Data Sheet

LTA4H

Human, recombinant, C-terminal FLAG-tag, and His-tag

Catalog #: 70001

Lot#: 140417 **Conc.:** 1.4 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 100 mM imidazole, and 20% glycerol.

Stability: >6 months at -80°C. Avoid repeated freeze/thaw.

References:

1. Snelgrove, R.J., *et al.*, *Science* 2010: **330(2000)**:90-94.
2. Stsiapanava, A., *et al.*, *PNAS* 2014: **111(11)**:4227-4232.

Description: Human leukotriene A-4 hydrolase, also known as LTA4H, LTA4, and LTA4 hydrolase (GenBank Accession No. NM_000895, a.a. 1-611(end) with C-terminal FLAG-tag and His-tag, MW = 71 kDa, expressed in an *E. coli* cell expression system.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

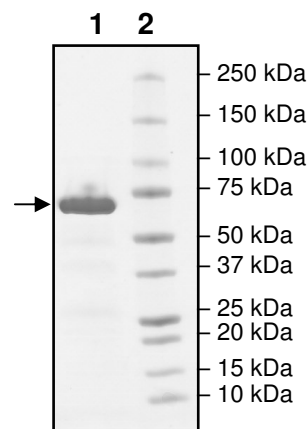
4-20% SDS-PAGE Coomassie staining

Lane 1:
2.8 µg LTA4H

Lane 2:
Protein Marker

Purity: ≥90%

MW: 71 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MPEIVDTCSLASPASVCRTKHLHLRCSVDFTRRTLGTAAALTVQSQEDNLRSLV
LDTKDLTIEKVINGQEVKYALGERQSYKGSPMEISLPIALSKNQEIVIEISFETSP
KSSALQWLTPEQTSKGHEPYLFSQCQAIHCRAILPCQDTPSVKLTYTAEVSVPK
ELVALMSAIRDGETPDPEDPSRKIYKFIQKVPIPCYLIALVVGALSRQIGPRTL
WSEKEQVEKSAYEFSETESMLKIAEDLGGPYVWGQYDLLVLPSPFPYGGMENP
CLTFVTPTLLAGDKSLSNVIAHEISHSWTGNLVTNKTWDHFWLNEGHTVYLERHI
CGRLFGEKFRHFNALGGWGELQNSVKTFGETHPFTKLVDLTDIDPDVAYSSV
PYEKGFA LLFYLEQLLGGPEIFLGFLKAYVEKFSYKSITDDWKDFLYSYFKDKV
DVLNQVDWNAWLYSPGLPPIKPNYDMTLTNACIALSQRWITAKEDDLNSFNATD
LKDLSSHQLNEFLAQLQRAPLPLGHIKRMQEVYNFNAINNSEIRFRWLRLCIQS
KWEDAIPALKMATEQGRMKFTRPLFKDLAAFDKSHDQAVRTYQEHKASMHP
VTAMLVGKDLKVD**DYKDDDDKHHHHH**

Black: LTA4H sequence

Red: FLAG-tag

Blue: His-tag

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com