
Mouse Lung

(Catalog No. X12010)

Concentration

500 µg protein in 200 µl SDS-PAGE Western blotting buffer

Size

500 µg

Storage Buffer

20% Glycerol, 250 mM Tris pH 6.8, 10% β-Mercaptoethanol

Storage

Extracts should be stored at -20°C and repeated freezing and thawing should be minimized

Extract Origin

Mouse lung tissue

Applications

Western blotting positive control; recommended use is 50 µg (20 µl) per lane; sample vial should be boiled once prior to use

Ordering Information:

Products	Size	Cat. No.
Mouse Lung	500 µg	X12010-1

This product is for research purposes only. Not intended for use in diagnostic procedures.