

---

## Mouse Cerebellum

(Catalog No. X12008)

---

### 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 cerebellum 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 Cerebellum	500 µg	X12008-1

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