

**Hemin (porcine)**

sterile 500x stock solution in 30% triethanolamine

Cat. No.	Amount
ML-108S	2 ml (for 1 L medium)
ML-108L	10 ml (for 5 L medium)
ML-108XL	20 ml (for 10 L medium)
ML-108XXL	5x 20 ml (for 50 L medium)

For research use only! Not intended for human or animal diagnostic or therapeutic uses.

Shipping: shipped on gel packs

Storage Conditions: store at 4 °C

Additional Storage Conditions: store dark

Shelf Life: 12 months

CAS#: 16009-13-5 (Hemin), 102-71-6 (Triethanolamine)

EC number: 240-140-1 (Hemin), 203-049-8 (Triethanolamine)

Description:

0,25% (w/v) solution of porcine Hemin in 30% Triethanolamine, sterile. For preparation of growth media for *Leishmania tarentolae*.

Preparation of *Leishmania* cultivation media:

Add to 1 l medium:

2.0 ml Hemin stock solution (500x) to a final concentration of 5 µg/ml.