

**L-Glutathion reduced - 0.16 M**

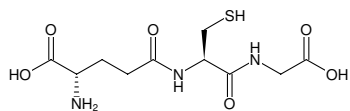
sterile filtered

 γ -L-Glutamyl-L-cysteinyl-glycine

Cat. No.	Amount
CSS-199	100 ml

Description:

Crystallization Stock Solution

**For *in vitro* use only!****Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Shelf Life:** 12 months**Molecular Formula:** $\text{H}_2\text{NCH}(\text{CO}_2\text{H})\text{CH}_2\text{CH}_2\text{CONHCH}(\text{CH}_2\text{SH})\text{CONHCH}_2\text{CO}_2\text{H}$ **Molecular Weight:** 307.32 g/mol**CAS#:** 70-18-8**EC number:** 200-725-4**Form:** liquid**Concentration:** 0.16 M