

---

## DTT Solution

(Catalog No. R-1104)

---

### Description

DTT, also known as Cleland's Reagent, is a particularly strong reducing agent, owing to its high conformational propensity to form a six member ring with an internal disulfide bond. It has a redox potential of -0.33 V @ pH 7. It is used for many protein research applications, including reducing polyacrylamide electrophoresis (SDS-PAGE). The 500mM stock solution can be directly used to achieve the amount of reducing agent required for your specific application.

### Molecular Weight (kDa)

154 kDa

### Formulation

10 mM phosphate buffered saline, pH 7.4

### Concentration/Volume

500 mM, 1 ml

### Storage Temp & Conditions

>6 months at 4°C.

### Application Notes

Used in biochemistry works to reduce disulfide bridges, protect biomolecules, in sample preparation, and to denature proteins before electrophoresis analysis (SDS-PAGE). 1-10 mM to maintain reduced proteins in solution, 50-100 mM for complete reduction for electrophoresis.

### Ordering Information:

Products	Size	Cat. No.
DTT Solution	500 mM, 1 ml	R-1104 -1

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