

**BIS-TRIS - 1 M**

any pH value

useful buffer range: pH 5.8 - 7.2

2-Bis (2-hydroxyethyl) amino-2-(hydroxymethyl)-1,3-propanediol

Cat. No.	Amount
CSS-503	100 ml

**Description:**

Crystallization Stock Solution, 1M BIS-TRIS adjusted with HCl, sterile filtered

**For in vitro use only!****Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Shelf Life:** 12 months**Molecular Formula:** C<sub>8</sub>H<sub>19</sub>NO<sub>5</sub>**Molecular Weight:** 209.24 g/mol**CAS#:** 6976-37-0**EC number:** 230-237-7**Form:** liquid**Concentration:** 1 M**pH:** arbitrary in 0.1 increments