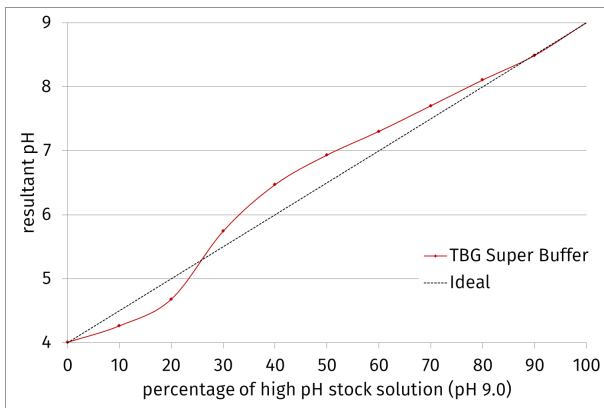




## TBG pH 9.0 - 1 M

Buffer System: Sodium Tartrate Dihydrate, BIS-TRIS and Glycylglycine  
sterile filtered

Cat. No.	Amount
CSS-385	100 ml



TBG Super Buffer: Mixing low (CSS-384) and high (CSS-385) pH stock solutions results in an almost linear pH function

**For *in vitro* use only!**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at 4 °C

**Shelf Life:** 12 months

**Molecular Formula:**

$\text{Na}_2\text{C}_4\text{H}_4\text{O}_6 \times 2 \text{H}_2\text{O}$

$\text{C}_8\text{H}_{19}\text{NO}_5$

$\text{C}_4\text{H}_8\text{N}_2\text{O}_3$

**Molecular Weight:** 230.08 g/mol; 209.24 g/mol; 132.12 g/mol

**CAS#:** 6106-24-7, 6976-37-0, 556-50-3

**Form:** liquid

**Concentration:** 1 M

**pH:** 9.0, adjusted with NaOH

**Description:**

Crystallization Stock Solution - TBG Buffer System: Sodium Tartrate Dihydrate, BIS-TRIS and Glycylglycine (3:2:2 Molar Ratio), pH 9.0 adjusted with NaOH

*Sodium Tartrate Dihydrate*

**Molecular Formula:**  $\text{Na}_2\text{C}_4\text{H}_4\text{O}_6 \times 2 \text{H}_2\text{O}$

**Molecular Weight:** 230.08

**CAS Number:** 6106-24-7

*BIS-TRIS*

**Molecular Formula:**  $\text{C}_8\text{H}_{19}\text{NO}_5$

**Molecular Weight:** 209.24

**CAS Number:** 6976-37-0

*Glycylglycine*

**Molecular Formula:**  $\text{C}_4\text{H}_8\text{N}_2\text{O}_3$

**Molecular Weight:** 132.12

**CAS Number:** 556-50-3

**Related Products:**

TBG pH 4.0, #CSS-384