

**PBS Buffer pH 7.6 - 10x conc.**

Phosphate buffered saline Buffer

Cat. No.	Amount
BU-115S-10	100 ml
BU-115L-10	1 l

**For *in vitro* use only!****Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Shelf Life:** 12 months**Form:** liquid**Concentration:** 10x**pH:** 7.6**Description:**

A 1 x concentration of PBS (Phosphate buffered saline) Buffer contains:

8.01 g/l NaCl

0.2 g/l KCl

1.78 g/l Na<sub>2</sub>HPO<sub>4</sub>·2H<sub>2</sub>O0.27 g/l KH<sub>2</sub>PO<sub>4</sub>

The Buffer is used for cell culture applications since it is isotonic and non-toxic to cells.

**Quality:**

- Molecular biology grade
- sterile
- DNase / RNase free