

Optiblot Buffer Recipes

Optiblot Reducing SDS Run Buffer (20X) – ab119195

- 0.6 M MOPS
- 1.2 M Tris
- 2% SDS
- 130 mM Sodium Bisulfite

MOPS (free acid)	62.80 g
Tris (free base)	72.60 g
SDS	10.0 g
Sodium Bisulfite	6.5 g
Ultra pure water to	500 ml (~385 g)

The pH should be between 8.2 and 8.3 @ 25 °C.

Optiblot SDS Run Buffer (20X) – ab119197

- 0.8 M Tricine
- 1.2 M Tris
- 2% SDS
- 50 mM Sodium Bisulfite

Tricine (free base)	71.70 g
Tris (free base)	72.60 g
SDS	10.0 g
Sodium Bisulfite	2.5 g
Ultra pure water to	500 ml (~385 g)

The pH should be between 8.2 and 8.3 @ 25 °C.

For Non-reduced samples (especially antibodies), omit the Sodium Bisulfite

For best results use low conductance ingredients to formulate the buffers.

Do NOT use acid or base to adjust pH of buffers.

Optiblot LDS Sample Buffer (4X) – ab119196

- 40% Glycerol
- 4% LDS, 0.8 M Triethanolamine-HCl pH 7.6
- 4% Ficoll-400
- 0.025% Phenol Red
- 0.025% Coomassie Brilliant Blue G250
- 2 mM EDTA-2Na

Glycerol	4.0 g
Triethanolamine	1.2 g
6N HCl	0.93 g
Lithium Dodecyl Sulfate	0.40 g
Ficoll 400	0.40 g
EDTA Di-Sodium	0.007 g
Brilliant Blue G250	0.0025 g
Phenol red	0.0025 g
Ultra pure water to	10 ml (~4.5 g)

The pH should be between 7.7 and 7.8 @ 25 °C.

For best results use low conductance ingredients to formulate the buffers.

Do NOT use acid or base to adjust pH of buffers.

Optiblot TGS Blot Stock Buffer (10X/20X) – ab119198

- 1.92 M Glycine
- 0.25 M Tris
- 1% SDS

Glycine	72.10g
Tris (free base)	15.20 g
SDS	5.0 g
Ultra pure water to	500 ml (~385 g)

The pH should be between 8.5 and 8.6 @ 25°C.

For best results use low conductance ingredients to formulate the buffers.

Do NOT use acid or base to adjust pH of buffers.

Instructions: Prior to use, dilute this blot buffer stock solution:

- 10X for use in semi-dry blotters
- 20X for use in Tank Blotters