
EpiQuik 6x Loading Dye Bromophenol Blue

(Catalog No. R12007)

Description

EpiQuik 6x Loading Dye Bromophenol Blue is used for loading samples on agarose or polyacrylamide gels.

Composition

- 10 mM TrisHCl (pH 7.6)
- 0.03% bromophenol blue
- 0.03% xylene cyanol FF
- 60% glycerol
- 60 mM EDTA

Quality Control

Functionally tested for DNA sample loading on agarose gel electrophoresis.

Storage Conditions

Store at 4°C up to 12 months.

Safety Warnings and Precautions

This product is designed for research purposes and in vitro use only. According to common laboratory safety practice, it is recommended to wear protective clothing, gloves and safety glasses. Please refer to www.epigentek.com for Material Safety Data Sheet of the product. Some applications this product is used in may require a license which is not provided by the purchase of this product. Users should obtain the license if required.

Ordering Information

Products	Size	Cat. No.
EpiQuik 6x Loading Dye Bromophenol Blue	6 x 1ml	R12007-1

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