



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

6X DNA Loading Buffers (Blue or Orange)

Loading buffers for gel electrophoresis containing density agents for loading DNA samples on gels and your choice of blue or orange tracking dyes.



Product attributes

Storage Conditions

Store at 2 to 8 °C or below

Product Description

6X DNA Loading Buffers contains density agents for loading DNA samples on agarose or polyacrylamide gels. Available with blue or orange tracking dyes.

- Density agents for easy sample loading
- Visible blue or orange dyes for tracking gel electrophoresis
- Compatible with agarose and polyacrylamide gels
- No fluorescent DNA stains

6X DNA Loading Buffer with Blue Tracking Dyes contains two blue tracking dyes that run at ~1.5 kb and ~200 bp in a 1% agarose gel. 6X DNA Loading Buffer (Orange) contains orange tracking dye that runs at ~50 bp in a 1% agarose gel. Unlike blue tracking dyes, orange dye does not shadow DNA bands, and is suitable for use with [DNAzure® Blue DNA Gel Stain](#).

These 6X DNA Loading Buffers do not contain fluorescent DNA dye. See our [GelRed® Prestain Plus 6X DNA Loading Dye](#) and [6X GelRed® Prestain Loading Buffers](#) for one-step loading and fluorescent staining. PCR products from [EvaGreen®](#) qPCR reactions also can be visualized on gels without additional fluorescent dye.

This datasheet was generated on June 8, 2026 at 07:37:29 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/6x-dna-loading-buffer-blue/>