



Glowing products for science

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

Go-Go™ Fast DNA Gel Running Buffer, 50X

A novel, low ionic strength running buffer that allows you to run DNA agarose gels at higher voltage to get results up to 3X faster than with TAE or TBE buffer.



Product attributes

Product Description

Go-Go™ Fast DNA Gel Running Buffer is a novel low ionic strength running buffer that allows you to run DNA agarose gels at higher voltage to get results up to 3X faster than with TAE or TBE buffer, streamlining routine DNA gel analysis. Go-Go™ Fast buffer also gives crisper band resolution compared to conventional electrophoresis buffers.

- Run your gel up to three times faster than with TAE or TBE buffer
- Crisp band resolution
- Excellent results with GelRed®, GelGreen®, and other popular dyes
- Convenient 50X concentrated stock solution

Run Your Gel 3 Times Faster with Go-Go™ Fast

Go-Go™ Fast DNA Gel Running Buffer gives excellent results with [GelRed® and GelGreen® DNA gel stains](#), as well as with [GelRed® Prestain Plus 6X DNA Loading Dye](#). The buffer can be used with other commonly used gel stains like SYBR® Safe and ethidium bromide. Go-Go™ Fast is supplied as a convenient 50X concentrated stock solution.

[View our full selection](#) of unique products for every step of your gel electrophoresis workflow, including DNA and RNA gel stains, DNA ladders, loading buffers, and more.

This datasheet was generated on January 9, 2026 at 11:52:15 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/go-go-fast-dna-gel-running-buffer-50x/>