

GelGreen™ Nucleic Acid Gel Stain FAQ

Question	Answer
Can GelGreen be used to stain ssDNA or RNA?	GelGreen can be used to stain ssDNA and RNA, but we recommend GelRed for this application because it is five times more sensitive for single stranded nucleic acids than GelGreen.
Is GelGreen compatible with downstream applications such as cloning, ligation and sequencing?	Yes. We recommend Qiagen or Zymoclean gel extraction kits or phenol-chloroform extraction to remove the dye from DNA.
Is GelGreen compatible with Southern or northern blotting?	GelGreen has not been validated in blotting applications.
Can I reuse a GelGreen precast gel after electrophoresis?	We do not recommend reusing GelGreen precast gels as signal decreases with subsequent electrophoresis.
How should I dispose of GelGreen?	GelGreen has passed the EPA regulated Title 22 test. Some facilities have approved the disposal of GelGreen directly down the drain. However, because regulations vary, please contact your safety office for local disposal guidelines. GelGreen can be adsorbed to activated carbon (also known as activated charcoal) for disposal as chemical waste.
What is the lower detection limit of GelGreen?	Some users have reported being able to detect bands containing less than 0.1 ng DNA. However, the limit of detection will depend on instrument capability and exposure settings.
Does GelGreen need to be used in the dark?	You can use the dye in room light, however we recommend storing the dye in the dark.
Is there a difference between 10,000X GelGreen in DMSO and water?	The GelGreen stock in water is a newer and improved product compared to the stock in DMSO. We recommend using GelGreen in water to avoid the potential hazards of handling DMSO, a solvent that can be absorbed through the skin. We continue to offer GelGreen in DMSO because some users do not wish to alter their established laboratory protocols.

Associated Products

GoldBio Catalog #	Product Name
A-201	Agarose LE (Molecular Biology Grade)
D010	1 kb DNA Ladder
D011	1 kb PLUS™ DNA Ladder
D001	100 bp DNA Ladder
P007	BLUestain™ Protein ladder, 11-245 kDa
P008	BLUestain™ 2 Protein ladder, 5-245 kDa
G-725	GelRed™ Nucleic Acid Stain Gel Stain, 10,000X in Water
E-670	EvaGreen® Dye, 20x in Water

GelGreen™ and its uses are covered by US patent numbers 7960498, 7803943, and 8232050. Materials from GoldBio are sold for research use only, and are not intended for food, drug, household, or cosmetic use.