



GenePath® Gel Kit

This GenePath Gel kit is composed of the following:

Item	Amount	Quantity
10 bottles, 100 ml 20x Running Buffer	100 ml	10
10 bottles, 100 ml 1% Agarose	100 ml	10
1% Low Melt Agarose and dropper	5 ml	1
1 mg/ml Ethidium Bromide and dropper	5 ml	1

Please refer to the GenePath reagent manual for more details.

1x Running Buffer

Pour 1 bottle of 20x running buffer into a 2 L graduated cylinder. QS to 2 L with distilled water.

Note: A precipitate may form upon prolonged storage. To redissolve, loosen bottle cap and autoclave for 15 minutes, or heat in a microwave until precipitate disappears.

1% Agarose

1. Melt a 100 ml bottle of agarose in a microwave. This will take 1 to 2 minutes. Check the agarose every 30 seconds to make sure the agarose does not boil over.
2. Pour the agarose into the GenePath casting stands. Let the agarose solidify for 30 to 60 minutes at room temperature before use.

1% Low Melt Agarose

1. Melt the low melt agarose in a microwave. This will take about 1 minute. Check the agarose every 15 seconds to make sure the agarose does not boil over.
2. This agarose is used to fasten the sample plug in the well.

1 mg/ml Ethidium Bromide

Ethidium bromide* is used to stain DNA fragments. Add 5 drops to approximately 300 ml distilled water. Always wear gloves when using ethidium bromide.

* Please refer to the MSDS.