

GoTaq® Flexi DNA Polymerase: Robust Performance with Magnesium Optimization

Article

Comments & Ratings

ABSTRACT

GoTaq® Flexi DNA Polymerase is a new member of the GoTaq® family of products. The two GoTaq® Flexi Buffers offer the flexibility of magnesium optimization in PCR. Green GoTaq® Flexi Buffer allows loading of amplified DNA samples directly onto agarose gels, eliminating the need to add sample loading buffer to PCR samples before electrophoresis. Colorless GoTaq® Flexi Buffer is provided for use when PCR samples are not loaded onto a gel or when the dyes in the Green GoTaq® Flexi Buffer would interfere with downstream applications.

Promega Notes 89, 17-20.

Kimberly Knoche, Becky Godat, Mark Koeff and Susan Fly Promega Corporation Publication Date: 2005



Download Article

(190 KB)



It appears that you have Javascript disabled. Our website requires Javascript to function correctly. For the best browsing experience, please enable Javascript.

1