

INCREASED STABILITY OF LONGLIFE™ ZYMOLYASE®

Enzymes regularly used in laboratory applications often require preparation of fresh solution before each use. Making fresh enzyme solution for each application is time consuming and wasteful. A wide variety of enzyme preparations in a ready-to-use format are offered, including LongLife™ Zymolyase®.

METHOD

LongLife™ Zymolyase® enzyme preparation was tested over a period of 4 weeks at 37°C and compared with LongLife™ Zymolyase® enzyme preparations stored at -20°C. LongLife™ Zymolyase® (1.5units/μl) was tested with freshly grown yeast suspension by monitoring the decrease in absorbance of the suspension.

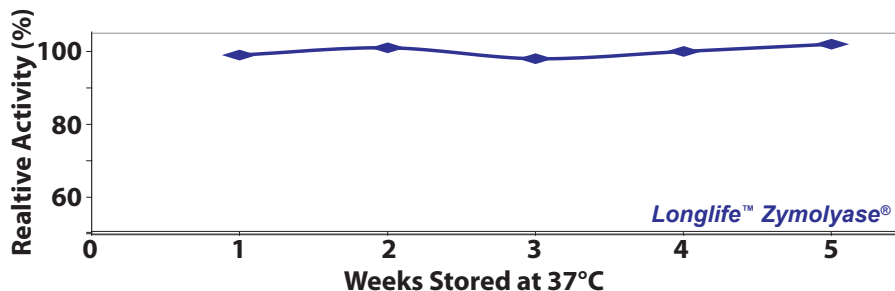


Figure 1: LongLife™ Zymolyase® is highly stable. LongLife™ Zymolyase® enzyme preparation was tested over a period of 4 weeks at 37°C and compared with LongLife™ Zymolyase® enzyme preparations stored at -20°C. LongLife™ Zymolyase® (1.5units/μl) was tested with freshly grown yeast suspension by monitoring the decrease in absorbance of the suspension. No measurable loss of activity was noticed.

CONCLUSION

No measurable loss of activity was noticed.

OTHER ENZYMES AVAILABLE

- LongLife™ Zymolyase® for the digestion of yeast and fungal cell walls.
- LongLife™ Lysozyme for the digestion of bacterial cell walls.
- LongLife™ PE LB Lysozyme for the digestion of bacterial cell walls and fully compatible with the PE LB buffer system. Reduces viscosity build-up due to presence of nucleases.
- LongLife™ Proteinase K for the digestion of proteins in nucleic acid preparations.
- LongLife™ Nuclease for the removal of nucleic acids.
- LongLife™ RNase for the digestion of RNA.
- LongLife™ DNase for the digestion of DNA.

REFERENCES

1. Gray, P. et al (2006) Mol Cell Prot. 6: 514
2. Saribas, A. et al (2004) Glycobiology 14: 1217



think proteins! think G-Biosciences!



G-Biosciences
St Louis, MO. 63132. USA
Phone: 1-800-628-7730
Fax: 1-314-991-1504
www.GBiosciences.com