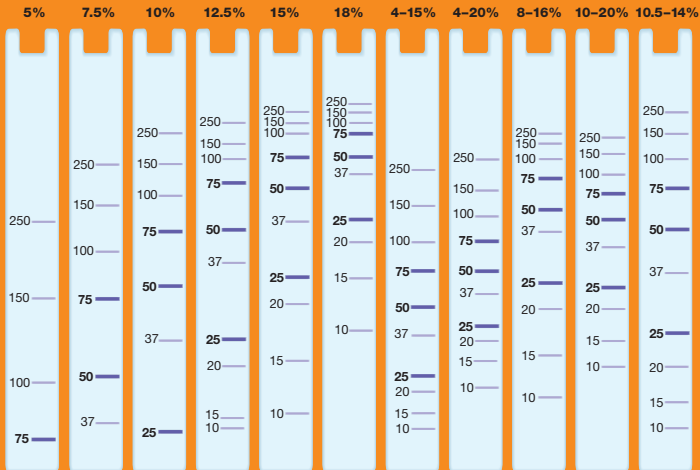


Tris-HCl Gel Protein Migration Chart

BIO-RAD

Precision Plus Protein™
Standards
Run on
Criterion™
Tris-HCl Gels

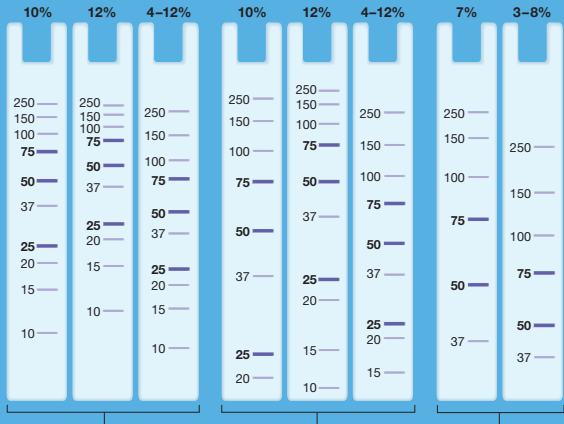


Tris-HCl gels with Tris/glycine/SDS running buffer

Neutral pH Gel Protein Migration Chart

BIO-RAD

Precision Plus Protein™ Standards
Run on Criterion™ XT Gels



Criterion XT Bis-Tris gels with XT MES running buffer: ideal for small proteins

Criterion XT Bis-Tris gels with XT MOPS running buffer: ideal for mid-size proteins

Criterion XT Tris-acetate gels with XT Tricine running buffer: ideal for large proteins

For ordering information, visit us on the Web at discover.bio-rad.com

NOTICE TO PURCHASER: LIMITED LICENSE

*This product is sold under license from Life Technologies Corporation, Carlsbad, CA for use only by the buyer of the product. The buyer is not authorized to sell or resell this product or its components.