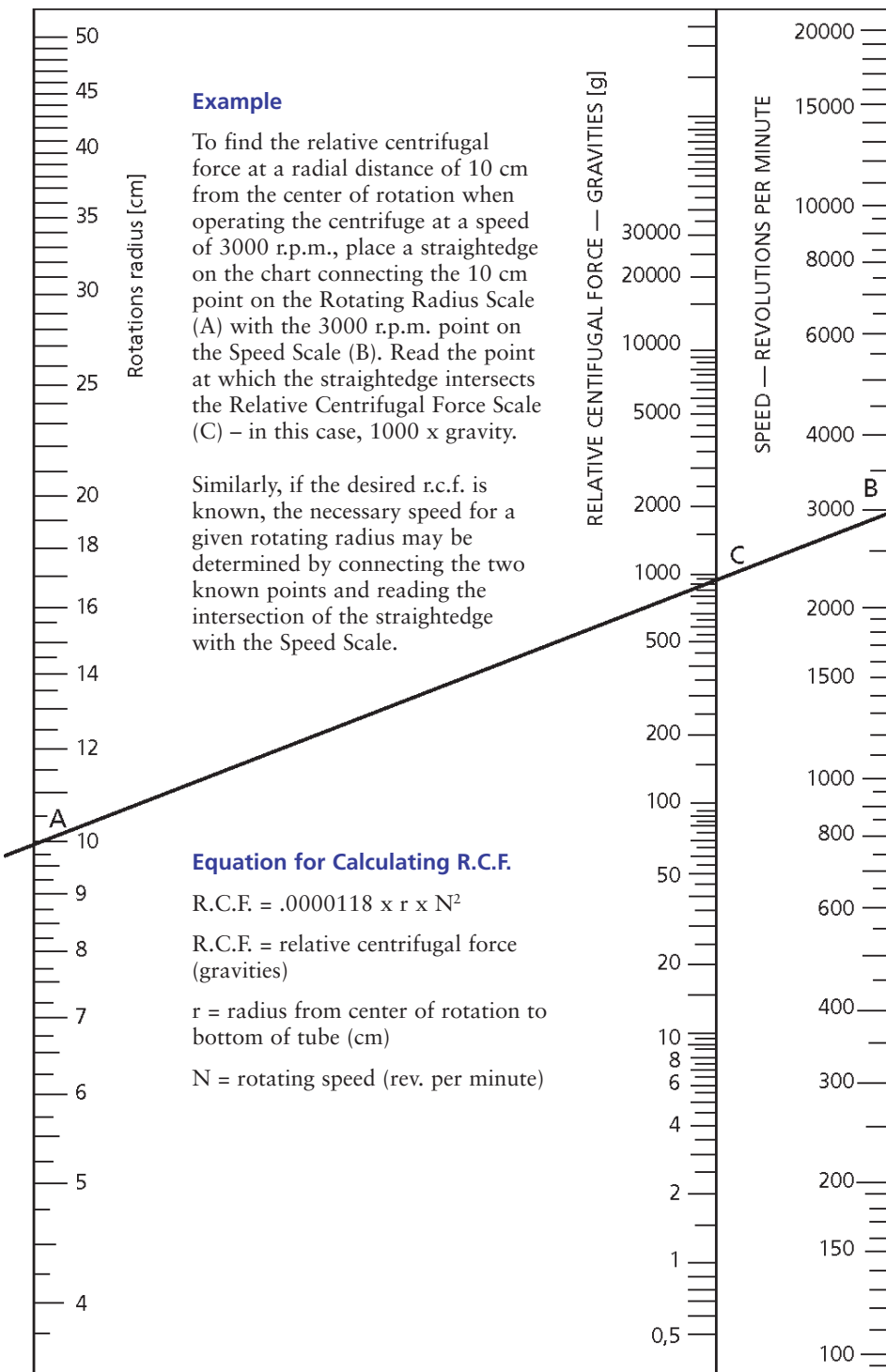


# Nomograph for RCF Determination



**BD Biosciences**

[bdbiosciences.com](http://bdbiosciences.com)

United States	Asia Pacific	Canada	Europe	Japan	Latin America/Caribbean
877.232.8995	65.6861.0633	888.259.0187	32.53.720.211	0120.8555.90	55.11.5185.9995

*For Research Use Only. Not for use in diagnostic or therapeutic procedures. Not for resale.*  
 BD, BD Logo and all other trademarks are the property of Becton, Dickinson and Company. ©2006 BD

