

Group 1: Red Blood Cells Products

Product No.	Product Name	Descriptions
816 056 100	Erytypecell A ₁ & B	A ₁ & B cells for ABO testing (2 x 10 mL) (<i>TANGO[®] optimo System use only</i>)
816 057 100	Biotestcell [®] A ₁ & B	A ₁ & B cells for ABO testing (2 x 10 mL)
816 047 100	Biotestcell [®] A ₂	A ₂ cells for ABO testing (1 x 10 mL)
816 014 100	Biotestcell [®] 1 & 2	Two cell antibody screen (2 x 10 mL)
816 085 100	Biotestcell [®] 3	Three cell antibody screen (3 x 10 mL)
816 022 100	Biotestcell [®] -I 11 Plus	Eleven cell antibody identification panel
816 065 100	Biotestcell [®] Pool	Pooled screening cells for donor antibody screening (1 x 10 mL)

2011 Schedule	Order by Date	Ship by Date	Expiration Date
	Jan-05-11	Feb-03-11	Mar-08-11
	Feb-02-11	Mar-03-11	Apr-05-11
	Mar-02-11	Mar-31-11	May-03-11
	Mar-30-11	Apr-28-11	May-31-11
	Apr-27-11	May-26-11	Jun-28-11
	May-25-11	Jun-23-11	Jul-26-11
	Jun-22-11	Jul-21-11	Aug-23-11
	Jul-20-11	Aug-18-11	Sep-20-11
	Aug-17-11	Sep-15-11	Oct-18-11
	Sep-14-11	Oct-13-11	Nov-15-11
	Oct-12-11	Nov-10-11	Dec-13-11
	Nov-09-11	Dec-08-11	Jan-10-12
Dec-07-11	Jan-05-12	Feb-07-12	

2012 Schedule	Order by Date	Ship by Date	Expiration Date
	Jan-04-12	Feb-02-12	Mar-06-12
	Feb-01-12	Mar-01-12	Apr-03-12
	Feb-29-12	Mar-29-12	May-01-12
	Mar-28-12	Apr-26-12	May-29-12
	Apr-25-12	May-24-12	Jun-26-12
	May-23-12	Jun-21-12	Jul-24-12
	Jun-20-12	Jul-19-12	Aug-21-12
	Jul-18-12	Aug-16-12	Sep-18-12
	Aug-15-12	Sep-13-12	Oct-16-12
	Sep-12-12	Oct-11-12	Nov-13-12
	Oct-10-12	Nov-08-12	Dec-11-12
	Nov-07-12	Dec-06-12	Jan-08-13
Dec-05-12	Jan-03-13	Feb-05-13	

2013 Schedule*	Order by Date	Ship by Date	Expiration Date
	Jan-02-13	Jan-31-13	Mar-05-13
	Jan-30-13	Feb-28-13	Apr-02-13
	Feb-27-13	Mar-28-13	Apr-30-13
	Mar-27-13	Apr-25-13	May-28-13
	Apr-24-13	May-23-13	Jun-25-13
	May-22-13	Jun-20-13	Jul-23-13
	Jun-19-13	Jul-18-13	Aug-20-13
	Jul-17-13	Aug-15-13	Sep-17-13
	Aug-14-13	Sep-12-13	Oct-15-13
	Sep-11-13	Oct-10-13	Nov-12-13
	Oct-09-13	Nov-07-13	Dec-10-13
	Nov-06-13	Dec-05-13	Jan-07-14
Dec-04-13	Jan-02-14	Feb-04-14	

2014 Schedule*	Order by Date	Ship by Date	Expiration Date
	Jan-01-14	Jan-30-14	Mar-04-14
	Jan-29-14	Feb-27-14	Apr-01-14
	Feb-26-14	Mar-27-14	Apr-29-14
	Mar-26-14	Apr-24-14	May-27-14
	Apr-23-14	May-22-14	Jun-24-14
	May-21-14	Jun-19-14	Jul-22-14
	Jun-18-14	Jul-17-14	Aug-19-14
	Jul-16-14	Aug-14-14	Sep-16-14
	Aug-13-14	Sep-11-14	Oct-14-14
	Sep-10-14	Oct-09-14	Nov-11-14
	Oct-08-14	Nov-06-14	Dec-09-14
	Nov-05-14	Dec-04-14	Jan-06-15
Dec-03-14	Jan-01-15	Feb-03-15	

*Estimated date.

Bio-Rad Customer Service
 1-800-2BIO-RAD (1-800-224-6723) | Fax: 510-741-6373 | diagcs@bio-rad.com



Group 2: Red Blood Cells Products

Product Codes	Product Name	Descriptions
816 020 100	Biotestcell®-I 8	Eight cell antibody identification panel
816 021 100	Biotestcell®-I 11	Eleven cell antibody identification panel
816 030 100	Coombscell-E	IgG coated red cells

2011 Schedule	Order by Date	Ship by Date	Expiration Date
	Jan-14-11	Feb-17-11	Mar-22-11
	Feb-11-11	Mar-17-11	Apr-19-11
	Mar-11-11	Apr-14-11	May-17-11
	Apr-08-11	May-12-11	Jun-14-11
	May-06-11	Jun-09-11	Jul-12-11
	Jun-03-11	Jul-07-11	Aug-09-11
	Jul-01-11	Aug-04-11	Sep-06-11
	Jul-29-11	Sep-01-11	Oct-04-11
	Aug-26-11	Sep-29-11	Nov-01-11
	Sep-23-11	Oct-27-11	Nov-29-11
	Oct-21-11	Nov-22-11	Dec-27-11
	Nov-18-11	Dec-22-11	Jan-24-12
Dec-16-11	Jan-19-12	Feb-21-12	

2012 Schedule	Order by Date	Ship by Date	Expiration Date
	Jan-13-12	Feb-16-12	Mar-20-12
	Feb-10-12	Mar-15-12	Apr-17-12
	Mar-09-12	Apr-12-12	May-15-12
	Apr-06-12	May-10-12	Jun-12-12
	May-04-12	Jun-07-12	Jul-10-12
	Jun-01-12	Jul-05-12	Aug-07-12
	Jun-29-12	Aug-02-12	Sep-04-12
	Jul-27-12	Aug-30-12	Oct-02-12
	Aug-24-12	Sep-27-12	Oct-30-12
	Sep-21-12	Oct-25-12	Nov-27-12
	Oct-19-12	Nov-20-12	Dec-25-12
	Nov-16-12	Dec-20-12	Jan-29-13
Dec-14-12	Jan-24-13	Feb-26-13	

2013 Schedule*	Order by Date	Ship by Date	Expiration Date
	Jan-11-13	Feb-14-13	Mar-19-13
	Feb-08-13	Mar-14-13	Apr-16-13
	Mar-08-13	Apr-11-13	May-14-13
	Apr-05-13	May-09-13	Jun-11-13
	May-03-13	Jun-06-13	Jul-09-13
	May-31-13	Jul-02-13	Aug-06-13
	Jun-28-13	Aug-01-13	Sep-03-13
	Jul-26-13	Aug-29-13	Oct-01-13
	Aug-23-13	Sep-26-13	Oct-29-13
	Sep-20-13	Oct-24-13	Nov-26-13
	Oct-18-13	Nov-19-13	Dec-24-13
	Nov-15-13	Dec-19-13	Jan-21-14
Dec-13-13	Jan-16-14	Feb-18-14	

2014 Schedule*	Order by Date	Ship by Date	Expiration Date
	Jan-10-14	Feb-13-14	Mar-18-14
	Feb-07-14	Mar-13-14	Apr-15-14
	Mar-07-14	Apr-10-14	May-13-14
	Apr-04-14	May-08-14	Jun-10-14
	May-02-14	Jun-05-14	Jul-08-14
	May-30-14	Jul-01-14	Aug-05-14
	Jun-27-14	Jul-31-14	Sep-02-14
	Jul-25-14	Aug-28-14	Sep-30-14
	Aug-22-14	Sep-25-14	Oct-28-14
	Sep-19-14	Oct-23-14	Nov-25-14
	Oct-17-14	Nov-18-14	Dec-23-14
	Nov-14-14	Dec-18-14	Jan-20-15
Dec-12-14	Jan-15-15	Feb-17-15	

*Estimated date.

© 2011 Bio-Rad Laboratories, Inc. 09/11 DG11-0652 U-116

Bio-Rad Customer Service
 1-800-2BIO-RAD (1-800-224-6723) | Fax: 510-741-6373 | diagcs@bio-rad.com

