

Reagent Red Blood Cell Products 2011 Shipping Schedule – Group 1

Product Codes	Product Name	Descriptions	Uses	
			Traditional Testing	TANGO optimo
816 056 100	Erytypecell A ₁ & B	A ₁ & B cells for ABO testing (2 x 10 mL)		X
816 057 100	Biotestcell® A ₁ & B	A ₁ & B cells for ABO testing (2 x 10 mL)	X	
816 047 100	Biotestcell® A ₂	A ₂ cells for ABO testing (1 x 10 mL)	X	
816 014 100	Biotestcell® 1 & 2	Two cell antibody screen (2 x 10 mL)	X	X
816 085 100	Biotestcell® 3	Three cell antibody screen (3 x 10 mL)	X	X
816 065 100	Biotestcell® Pool	Pooled screening cells for donor antibody screening (1 x 10 mL)		X

Note: Using the below shipping dates, reagent red cell products will have approximately five weeks shelf life upon delivery. Non-red blood cell reagents, which are part of these standing orders, may ship separately for delivery prior to the red cell ship by date.

Order by Date	Ship by Date	Expiration Date
January 5	February 3	March 8
February 2	March 3	April 5
March 2	March 31	May 3
March 30	April 28	May 31
April 27	May 26	June 28
May 25	June 23	July 26
June 22	July 21	August 23
July 20	August 18	September 20
August 17	September 15	October 18
September 14	October 13	November 15
October 12	November 10	December 13
November 9	December 8	January 10, 2012
December 7	January 5, 2012	February 7, 2012

See reverse side for **Group 2 Schedule** >

Bio-Rad Customer Service

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



Reagent Red Blood Cell Products

2011 Shipping Schedule – Group 2

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

Note: Using the below shipping dates, reagent red cell products will have approximately five weeks shelf life upon delivery. Non-red blood cell reagents, which are part of these standing orders, may ship separately for delivery prior to the red cell ship by date.

Order by Date	Ship by Date	Expiration Date
January 14	February 17	March 22
February 11	March 17	April 19
March 11	April 14	May 17
April 8	May 12	June 14
May 6	June 9	July 12
June 3	July 7	August 9
July 1	August 4	September 6
July 29	September 1	October 4
August 26	September 29	November 1
September 23	October 27	November 29
October 21	November 22	December 27
November 18	December 22	January 24, 2012
December 16	January 19, 2012	February 21, 2012

See reverse side for **Group 1 Schedule** >

Bio-Rad Customer Service

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

