



# Red Blood Cells Products – Group 1

## 2015 Schedule

Product No.	Product Name	Descriptions
816 056 100	Erytypecell A <sub>1</sub> & B	A <sub>1</sub> & B cells for ABO testing (2 x 10 mL) ( <i>TANGO® system use only</i> )
816 057 100	Biotestcell® A <sub>1</sub> & B	A <sub>1</sub> & B cells for ABO testing (2 x 10 mL)
816 047 100	Biotestcell® A <sub>2</sub>	A <sub>2</sub> 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)

Order by Date	Ship by Date	Expiration Date
Dec-10-2014	Jan-8-2015	Feb 9-2015
Jan-7-2015*	Feb 5-2015	Mar-16-2015
Feb-11-2015	Mar-12-2015	Apr-20-2015
Mar-11-2015	Apr-9-2015	May-18-2015
Apr-8-2015	May-7-2015	Jun-15-2015
May-6-2015	Jun-4-2015	Jul-13-2015
Jun-3-2015	Jul-2-2015	Aug-10-2015

Order by Date	Ship by Date	Expiration Date
Jul-1-2015	Jul-30-2015	Sep-7-2015
Jul-29-2015	Aug-27-2015	Oct-5-2015
Aug-26-2015	Sep-24-2015	Nov-2-2015
Sep-23-2015	Oct-22-2015	Nov-30-2015
Oct-21-2015**	Nov-19-2015	Dec-28-2015
Nov-11-2015	Dec-10-2015	Jan-18-2016
Dec-2-2015	Jan-7-2016	Feb-11-2016
Jan-6-2016	Feb-4-2016	Mar-15-2016

\*Five weeks until next shipment due to manufacturing schedule  
 \*\*Three weeks until next shipment due to manufacturing schedule

### Bio-Rad Customer Service

1-800-224-6723 | Fax: 510-741-6373 | [diagcs@bio-rad.com](mailto:diagcs@bio-rad.com)





# Red Blood Cells Products – Group 2

## 2015 Schedule

Product No.	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
* 816 500 100	Hemo-QC	Quality control for automation & manual testing

\*Schedule is for shipping information only. Expiration date will be provided with product.

Order by Date	Ship by Date	Expiration Date
Dec-24-2014	Jan-29-2015	Mar-9-2015
Jan-28-2015	Feb-26-2015	Apr-6-2015
Feb-25-2015	Mar-26-2015	May-4-2015
Mar-25-2015	Apr-23-2015	Jun-1-2015
Apr-22-2015	May-21-2015	Jun-29-2015
May-20-2015	Jun-18-2015	Jul-27-2015
Jun-17-2015	Jul-16-2015	Aug-24-2015

Order by Date	Ship by Date	Expiration Date
Jul-15-2015	Aug-13-2015	Sep-21-2015
Aug-12-2015	Sep-10-2015	Oct-19-2015
Sep-9-2015	Oct-8-2015	Nov-16-2015
Oct-7-2015*	Nov-5-2015	Dec-14-2015
Oct-28-2015	Nov-25-2015	Jan-4-2016
Nov-25-2015	Dec-24-2015	Feb-1-2016
Dec-30-2015	Jan-28-2016	Mar-8-2016

\*Three weeks between shipments due to manufacturing schedule.

### Bio-Rad Customer Service

1-800-224-6723 | Fax: 510-741-6373 | [diagcs@bio-rad.com](mailto:diagcs@bio-rad.com)

