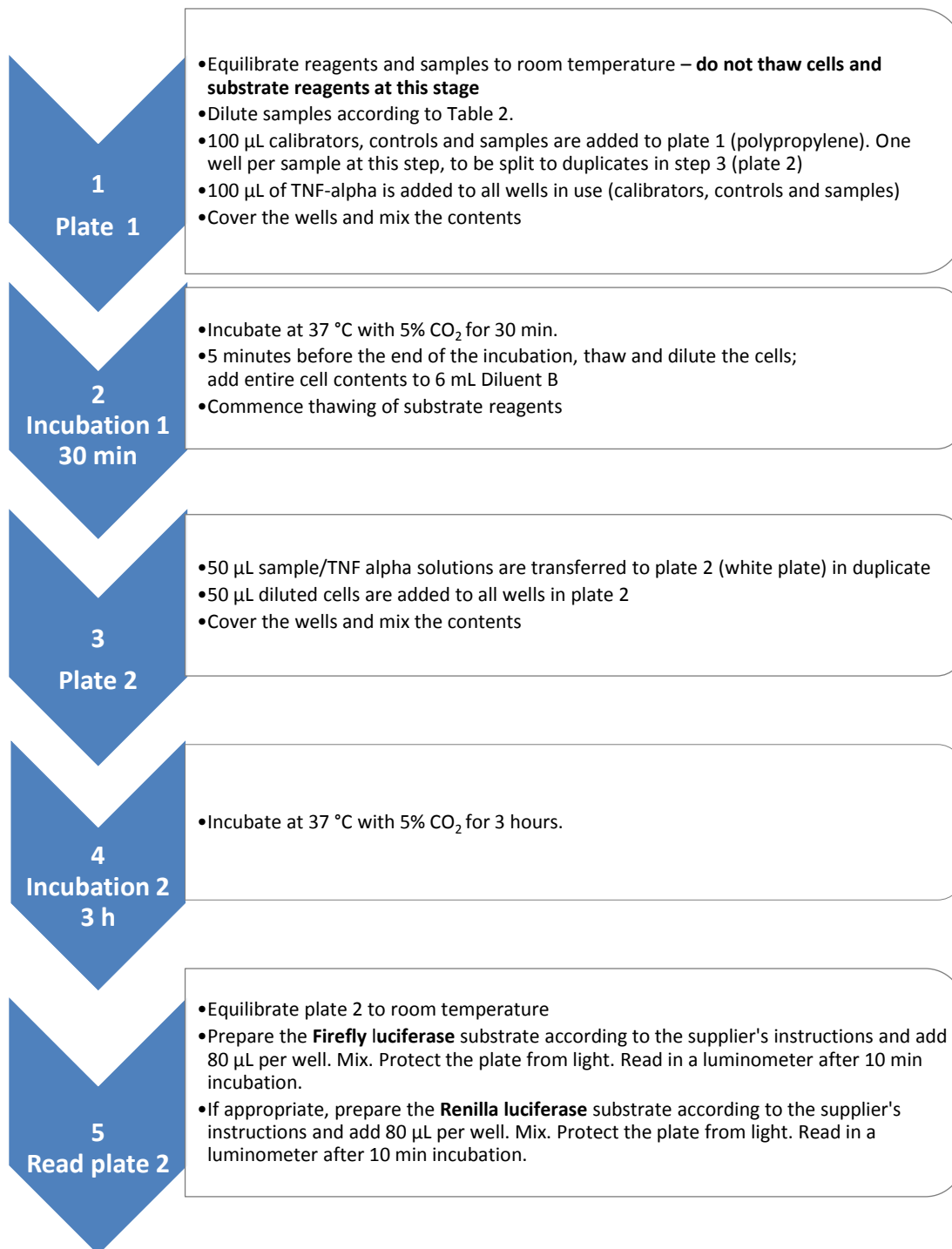




Quick Guide – quantification of TNF-alpha inhibitor biological activity



Troubleshooting, FAQ and Contacts

Please consult Euro Diagnostica's website www.eurodiagnostica.com

distributed in the US/Canada by:
Eagle Biosciences, Inc.
20A NW Blvd, Suite 112 Nashua, NH 03063
Phone: 617-419-2019 FAX: 617-419-1110
www.EagleBio.com info@EagleBio.com



EAGLE
BIOSCIENCES

Euro Diagnostica AB

Mail address:
P.O. Box 50117
SE - 202 11 Malmö
Sweden

Visiting address:
Lundavägen 151
Malmö
Sweden

T +46 40 53 76 00
F +46 40 43 22 88
E info@eurodiagnostica.com
W www.eurodiagnostica.com

Doc No: E-173-GB01, June 2016