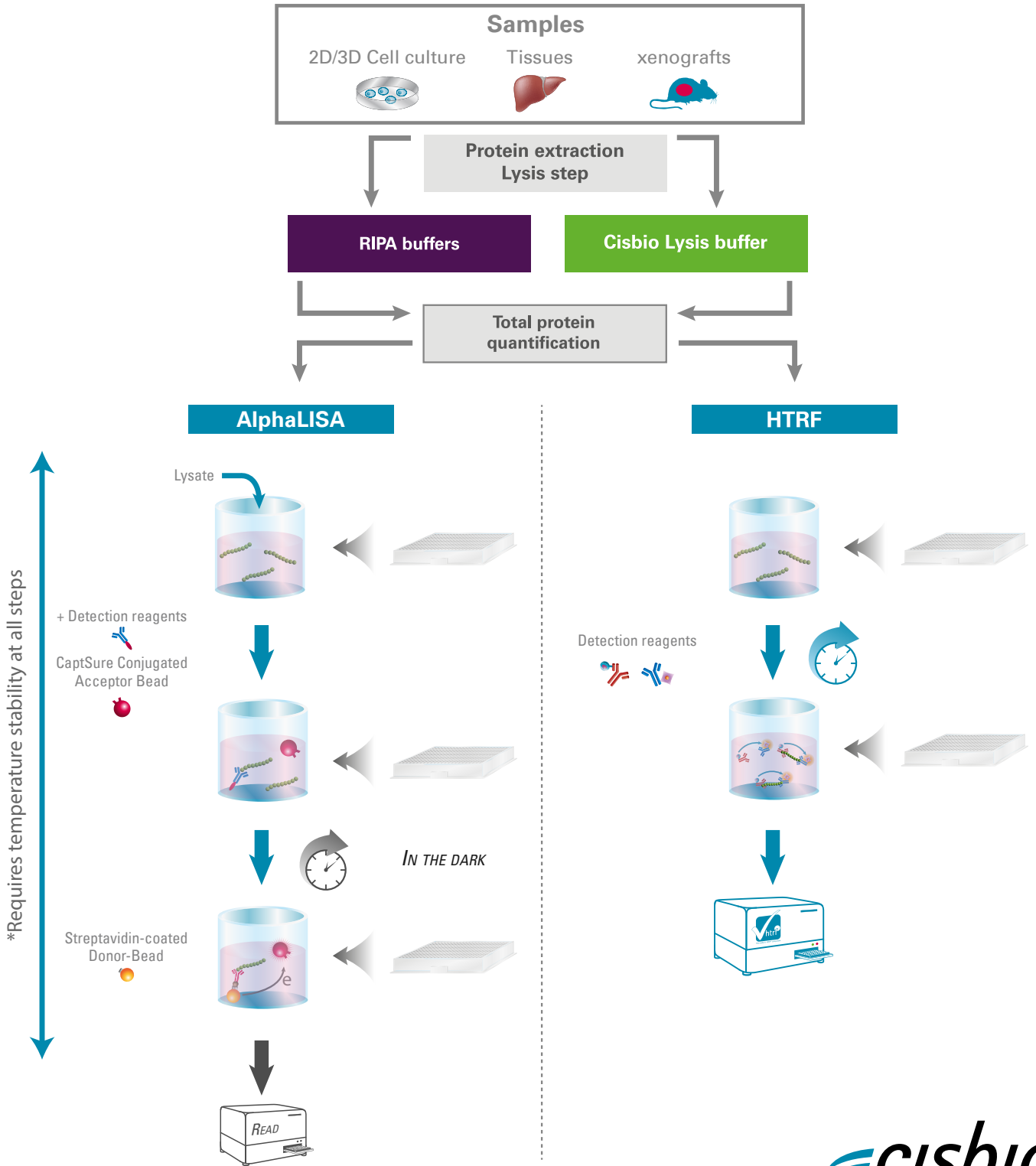




SAVE TIME AND MONEY

Switch to HTRF[®] assays



FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567 China +86 10 8080 9288

Visit www.cisbio.com to find a list of our regional distributors

