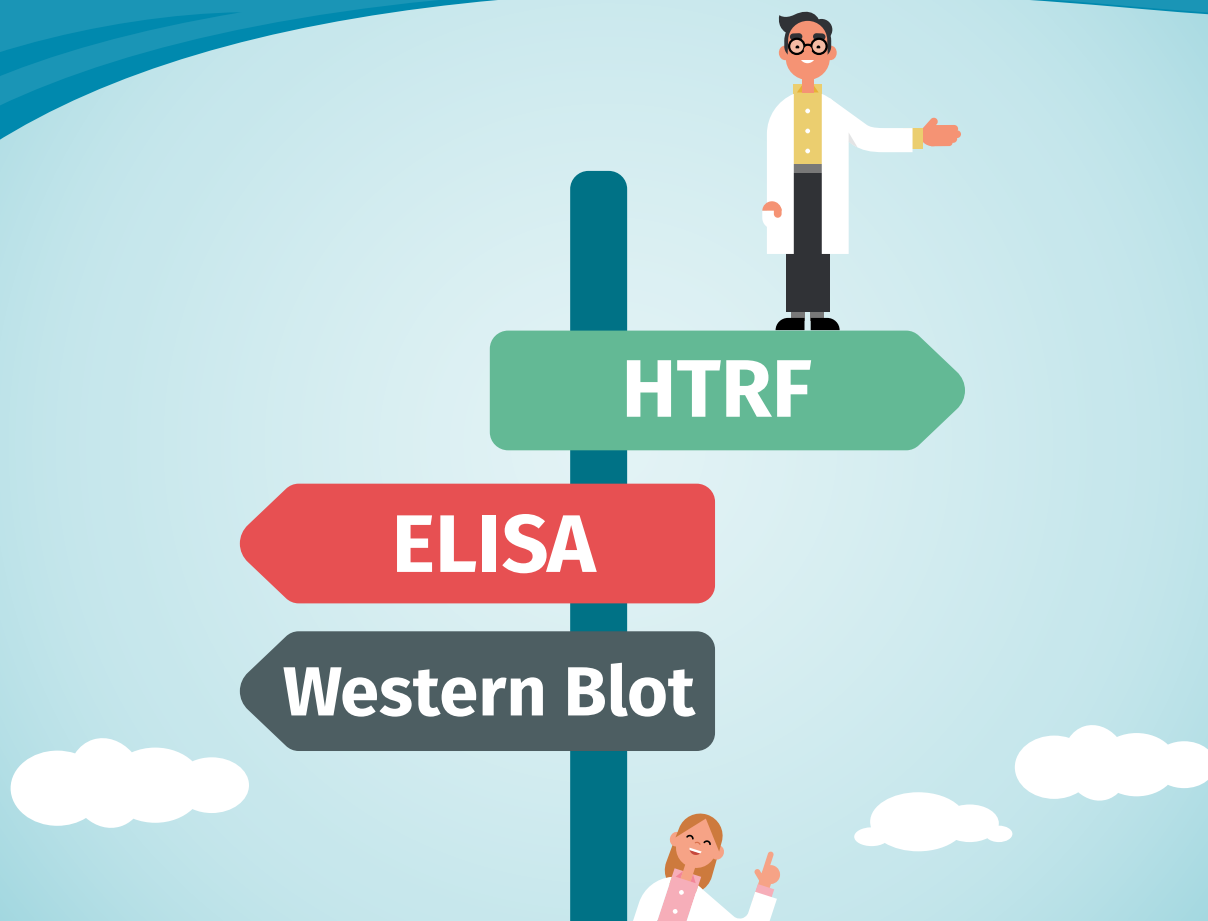


How **HTRF** compares to Western Blot and ELISA

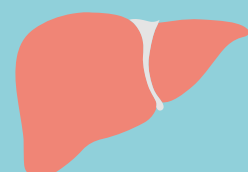


The path to excellent results does not have to be challenging.
Will you be convinced of the **ease** of **HTRF**?

LYSATE PREPARATION
from different type of samples:



2D/3D Cell culture

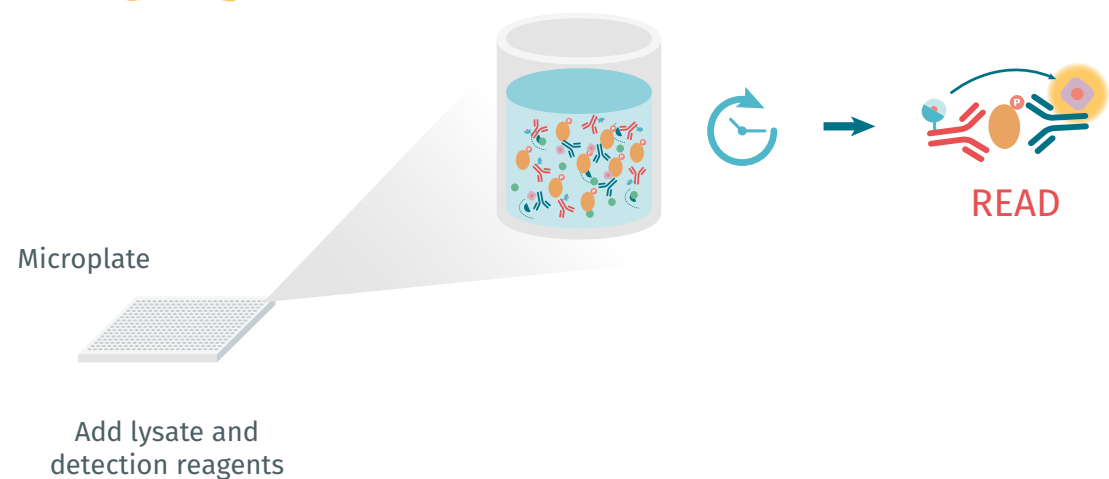


Tissues

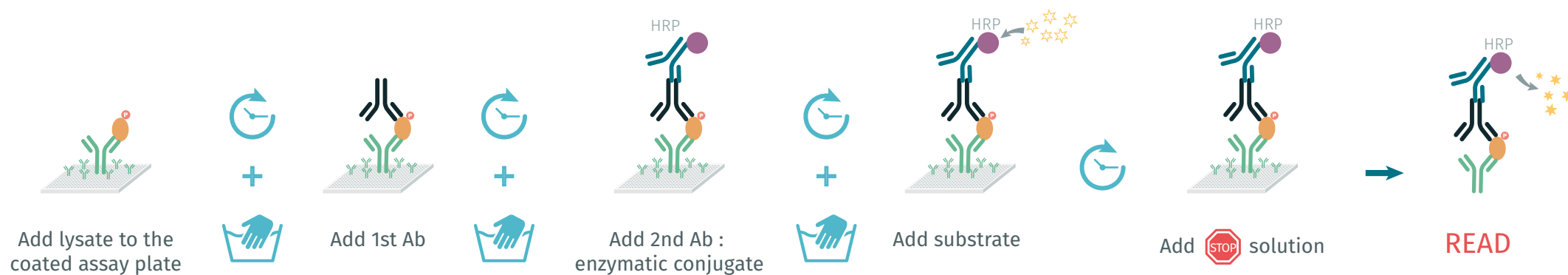


Xenografts

HTRF



ELISA

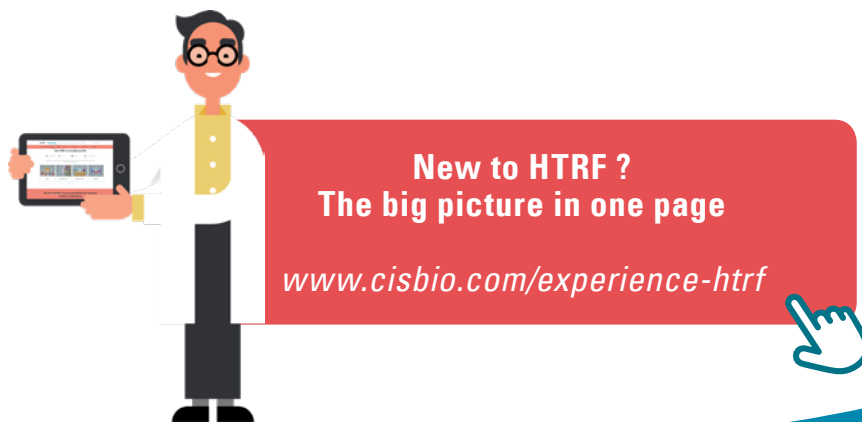
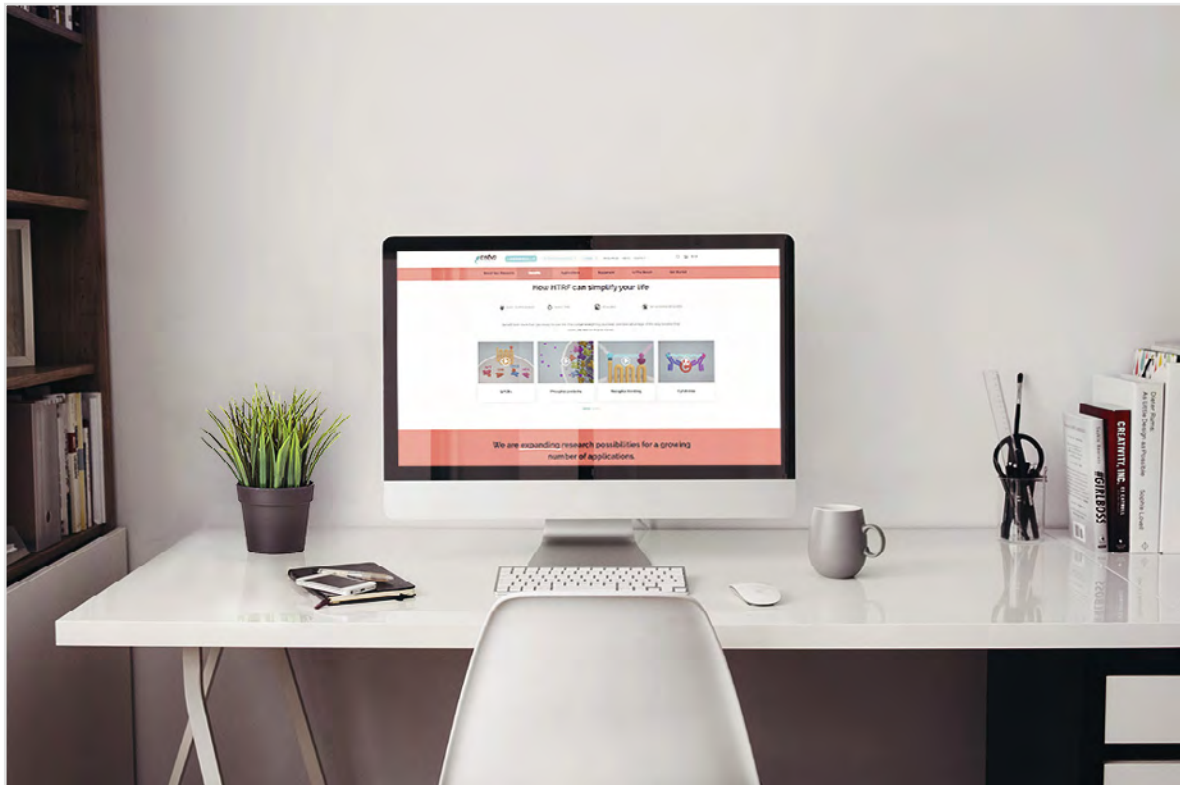


WESTERN BLOT



HTRF SIMPLIFY YOUR LIFE

Benefit from more than 300 ready-to-use kits that contain everything you need. Take advantage of this easy solution saving you both time and money.



FOR MORE INFORMATION

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

China +86 21 5018 9880 Japan +81-(0)43-306-8712

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

