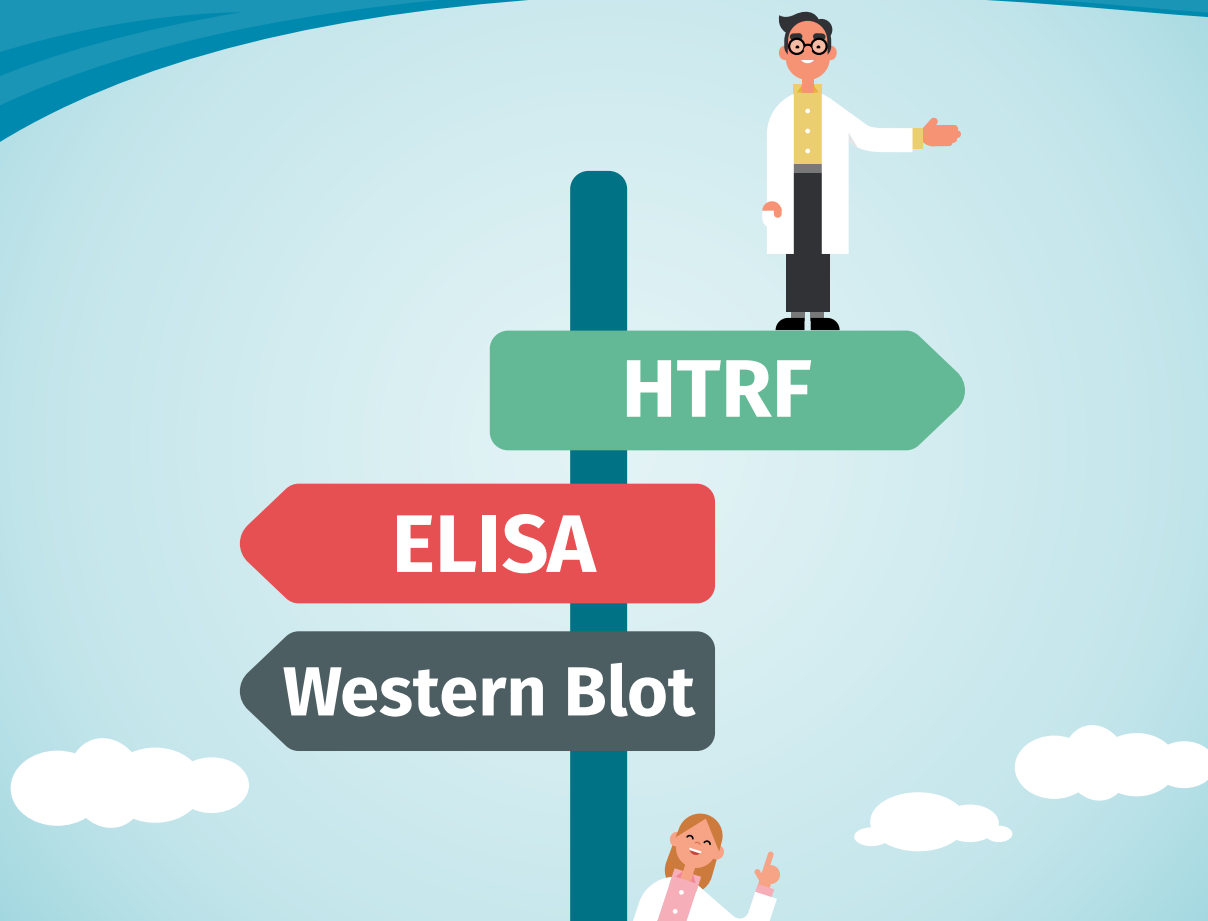


# How HTRF compares to Western Blot and ELISA



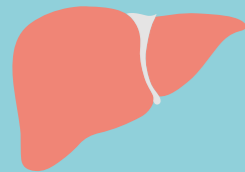
The path to excellent results does not have to be challenging.  
Will you be convinced of the **ease** of **HTRF**?

# LYSATE PREPARATION

from different type of samples:



2D/3D Cell culture



Tissues

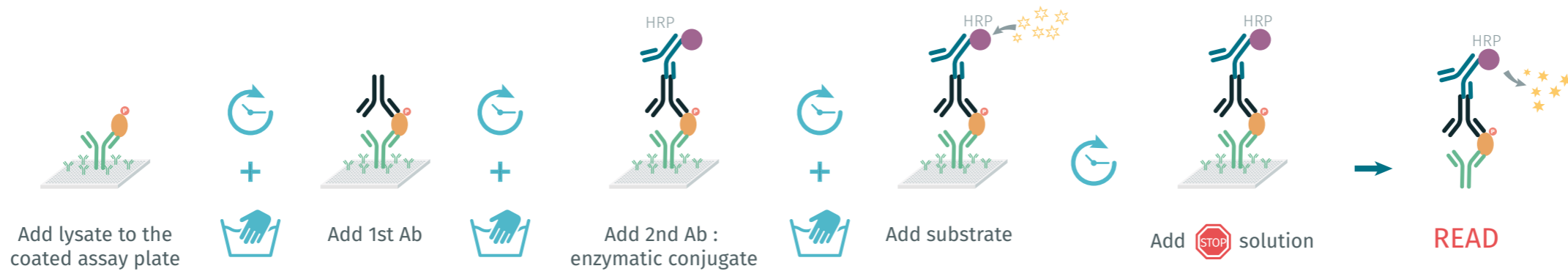


Xenografts

## 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.



**New to HTRF ?  
The big picture in one page**

[www.cisbio.com/experience-htrf](http://www.cisbio.com/experience-htrf)



## 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](http://www.cisbio.com) to find a list of our regional distributors



cisbio