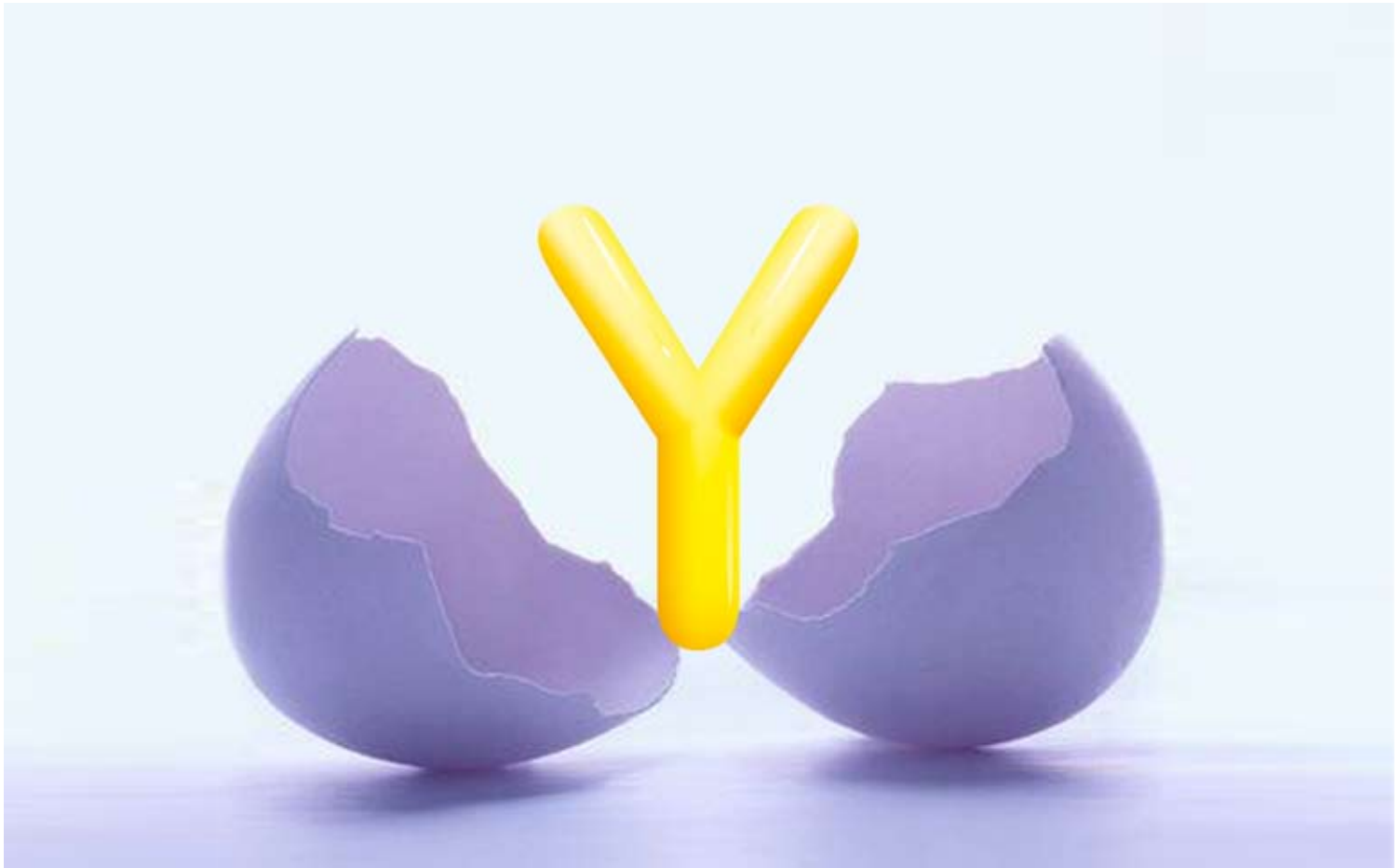


# AFFINITY PURIFIED CUSTOM CHICKEN ANTIBODIES (IgY)



WHY BLEED?

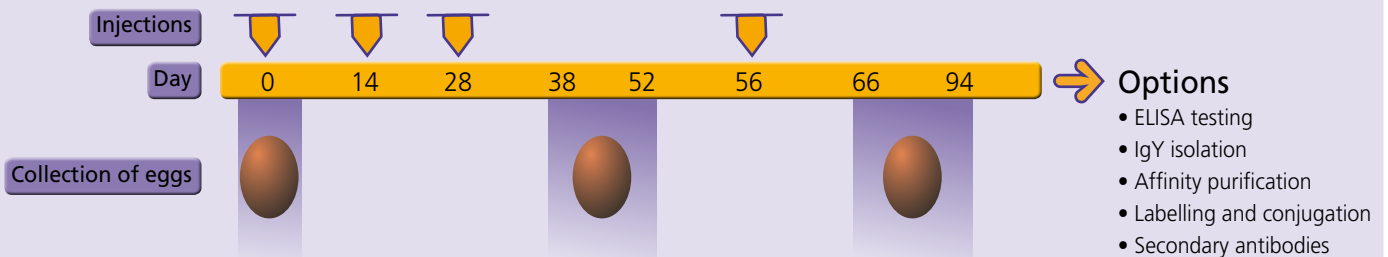
## WHY PRODUCING ANTIBODIES IN CHICKEN?

- **Improve your results**  
IgY Abs produce less background in your application.
- **Get rid of homology issues**  
Using a non-mammalian host for immunisation allows you to consider even highly conserved mammalian proteins as targets.
- **Obtain large Abs quantities**  
Only two eggs contain as much Antibody as in 100 ml of rabbit serum.
- **Get pure IgY antibodies...without trouble**  
Our team is used to perform IgY purification from egg yolk. Chose that option to save time and money!
- **Need secondary antibodies?**  
Anti-IgY secondary antibodies are available in our catalogue, with a rich variety of conjugates.

## FOR YOUR SPECIFIC NEEDS, EUROGENTEC OFFERS...

- > The production of antibodies in chicken
- > The isolation of IgY
- > The affinity purification of the target-specific antibody sub-population.

↓ Your protein of interest or peptide(s) designed by Eurogentec



## DID YOU KNOW THAT?

The antibody production in chicken is non-invasive, as the corresponding antibodies are not isolated from blood, but from eggs.

## How to start?

Contact us at [info@eurogentec.com](mailto:info@eurogentec.com) or visit our website [www.eurogentec.com](http://www.eurogentec.com).

We will design the best protocol and options corresponding to your needs.