



## Recombinant CloneDetect

An animal-free antibody reagent for detecting human IgG

### Key Features

- Free from raw materials of animal origin
- High specificity and affinity for all sub-classes of human IgG
- Specific to the Fc region
- Label-free, *in situ* detection of IgG secretion
- Each batch is subjected to rigorous safety assessment with full certification

Recombinant CloneDetect is a fluorescein-labeled antibody with high affinity and specificity for all sub-classes of human IgG. Recombinant CloneDetect is manufactured under animal-free methods in Chinese Hamster Ovary cells (CHO) using no raw materials of animal origin. Each manufactured batch of Recombinant CloneDetect is subjected to a rigorous safety assessment, with full certifications readily available upon request.

### Screen and select mammalian clones secreting human IgG

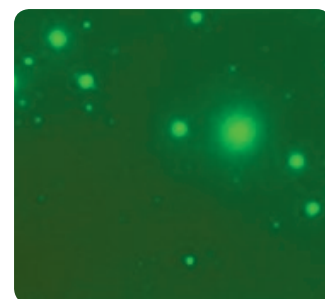
Recombinant CloneDetect is designed for use with ClonePix™ Systems and CloneMedia Semi-Solid Media when used to screen and select mammalian cell lines secreting human IgG. ClonePix Systems increase workflow productivity in over 100 biopharmaceutical laboratories around the world. Users can screen more clones in less time and select clones with optimal expression levels early in the workflow.

### Formation of discrete colonies

To prepare for screening, cells are grown in source plates containing semi-solid CloneMedia. This ensures formation of discrete colonies and facilitates plating out of large numbers of cells. Variants of the ready-to-use media are listed on the next page.

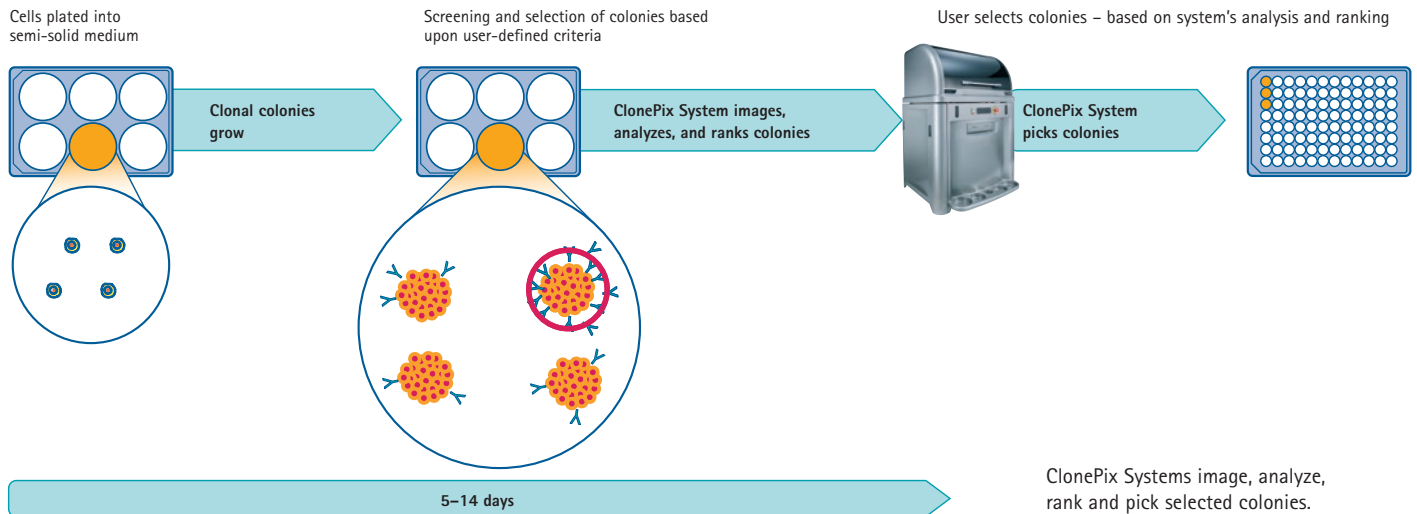
### Label-free, *in situ* detection of IgG secretion

Added to semi-solid CloneMedia at the moment of plating, Recombinant CloneDetect diffuses freely, creating immunoprecipitation complexes with target immunoglobulins as they are secreted from the colonies. These complexes generate localized fluorescent signals that, under optimized assay conditions, are proportional to the amount of IgG secreted by a specific clone. Utilizing a combination of imaging and robotics, ClonePix System screens for clonality and productivity simultaneously prior to collecting the highest producing clones in a rapid one-step process.



IgG secretion revealed: Fluorescent images of CHO-S cells taken on a ClonePix System 12 days after plating using Recombinant CloneDetect Media.

## Image, analyze, rank and pick secreting clones



### Supporting regulatory demands for minimized exposure to animal-derived components

Development and production of human IgG biotherapeutics must meet stringent regulatory demands, one of which is the ability to demonstrate minimal exposure to animal components throughout development and production. To support these regulatory requirements, Molecular Devices® offers a range of products free of raw materials of animal origin for use with the ClonePix Systems. All products are cell-culture validated, sterile and azide-free.

Visit [www.moleculardevices.com/genetix](http://www.moleculardevices.com/genetix) for further product information and the contact details of your local representative.

### Instrument compatibility



ClonePix 2 Systems

Recommended products for colony growth, human IgG detection and cell line expansion

Description	Part number
<b>Recombinant CloneDetect</b>	
Recombinant CloneDetect, anti-human IgG (Fc) Detection agent. Fluorescein labelled. Sterile, azide-free.	K8295
<b>CloneMedia and CloneMatrix</b>	
CloneMedia-CHO Semi-Solid Media for CHO-S cells	K8710
CloneMedia-CHO Semi-Solid Media for CHO-S cells (Bulk) • Contains (6) bottles of K8710	K8700
CloneMedia-CHO DHFR Semi-Solid Media for DG44, DUXB11 and CHO-S cells	K8712
CloneMedia-CHOK1SV Semi-Solid Media for CHOK1SV cells; glutamine free	K8725
CloneMedia-CHOK1SV Semi-Solid Media for CHOK1SV cells (Bulk); glutamine free • Contains (6) bottles of K8725	K8720
CloneMatrix Semi-Solid Media optimized for CHO cells, makes 100 mL final media volume*	K8530
CloneMatrix Semi-Solid Media (Bulk) optimized for CHO cells, makes 6 x 100 mL final media volume* • Contains (6) bottles of K8530	K8520
CloneMatrix Semi-Solid Media; makes 100 mL final media volume*	K8510
CloneMatrix Semi-Solid Media (Bulk); makes 6 x 100 mL final media volume* • Contains (6) of K8510	K8500

\* CloneMatrix is a 2.5x concentrate available for those who wish to make their own specific semi-solid media recipes.

#### Contact Us

Phone: +1-800-635-5577  
 Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
 Email: [info@moldev.com](mailto:info@moldev.com)  
 Check our website for a current listing of worldwide distributors.

#### Regional Offices

USA and Canada +1-800-635-5577  
 Brazil +55-11-3616-6607  
 China (Beijing) +86-10-6410-8669  
 China (Shanghai) +86-21-3372-1088  
 Germany 00800-665-32860

Japan (Osaka) +81-6-7174-8831  
 Japan (Tokyo) +81-3-6362-5260  
 South Korea +82-2-3471-9531  
 United Kingdom +44-118-944-8000

