

An Example of the Use of the iQue® High-Throughput Screening (HTS) by Cytometry Panel Builder

1

Select instrument by clicking on the dropdown box and then click next.

Cytometry Panel Builder

Step 1: Select Cytometer

Select Cytometer*

Type to search

- iQue 3 (BR)
- iQue 3 (VBR)
- iQue 3 (VYB)
- iQue 5 (VYBR)** Press enter to select
- iQue Screener PLUS (BR)
- iQue Screener PLUS (VBR)
- iQue Screener PLUS (VYB)

2

Enter panel requirements using the menu, including species reactivity, cell type, antibody vendor, or excluded fluorochromes/channels.

Cytometry Panel Builder

Step 2: Enter Panel Requirements

Species Reactivity *

Human

You can select species reactivity per each target from advanced settings later.

Sample Fixation

☒ No
Sample is not fixed or is fixed after staining.

☐ Yes
Sample is fixed before or during staining; Only compatible products/clones will be considered.

Sample Digestion (Collagenase)

☒ No
Sample is not prepared by digestion.

☐ Yes
Sample is prepared by digestion; Only compatible products/clones will be considered.

Cell Populations

Type to search

3

Select your markers and optionally assign fluorochromes.

Cytometry Panel Builder

Step 3: Enter Panel Details

Fill Panel Fields

Import Excel File Load Kits/Pre-defined Markers

1 Target / Marker* Fluorochrome / Special Dye Laser & Filter

Viability (Live/Dead, Live-Dead, L/D, ...) Live-Dead Fixable Violet Violet405 4...

2 Target / Marker* Expression Level Fluorochrome / Special Dye

CD3 (CD3e, CD3 epsilon, CD3 epsilon) High To Be Automatically Selected

3 Target / Marker* Expression Level Fluorochrome / Special Dye

CD4 High To Be Automatically Selected

4 Target / Marker* Expression Level Fluorochrome / Special Dye

CD8 High To Be Automatically Selected

4

Your panel will be generated and presented on this screen, here you will find a score of the suitability of the panel for your application, shown the spectral layout, and given more details about each antibody, including catalogue numbers for ordering.

Cytometry Panel Builder

Step 4: Get Optimized Panel Suggestion

Your Requested Panel EasyPanel Fully Computed Panel

iQue 5 (VYBR)

LASER	FILTER	TARGET	FLUOROCROME	PRODUCTS
Violet405	445/45	Viability (Live/Dead, Live-Dead, L/D, Live Dead)	* Live-Dead Fixable Violet	Select
Blue488	525/45	CD8	FITC	Select
Yellow561	586/20	CD25 (IL-2 Receptor alpha, IL-2 Receptor alpha, IL-2R alpha)	PE	Select
Yellow561	780/60	CD3 (CD3e, CD3 epsilon, CD3 epsilon, CD3e)	PE-Fix 810 (PE-FB10)	Select
Red637	725/40	CD4	Spark Red 718	Select

Edit Panel Manually

Share Export To Excel

Your EasyPanel Quality Score™

10/10 Details Spillover Matrix Balanced Thresholds Mode

EasyPanel Quality Score™ For Fully Computed Panel