

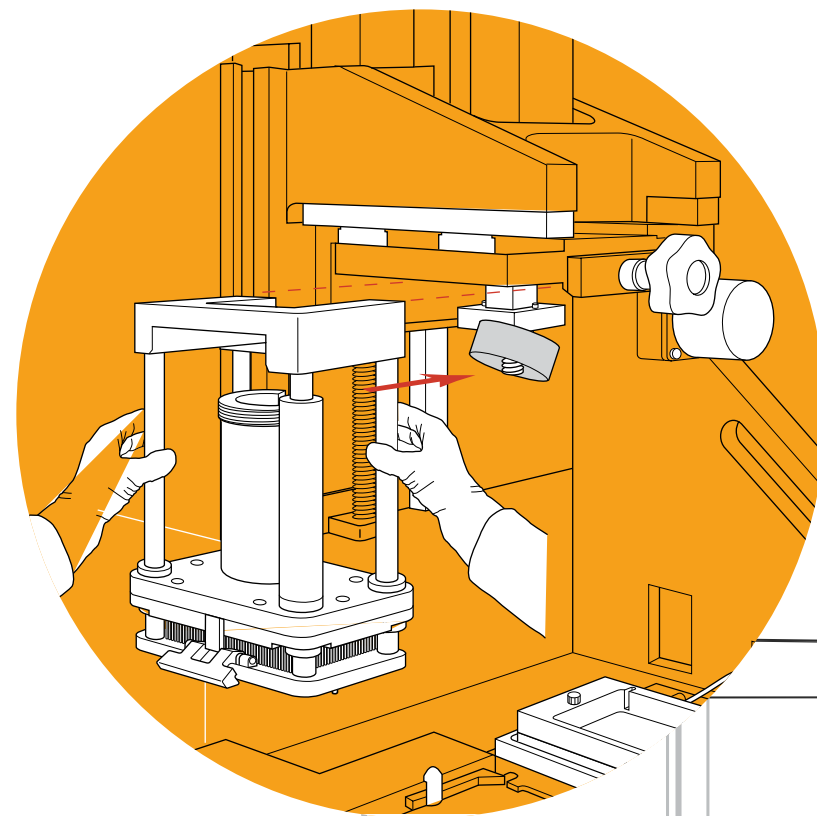
FLIPR Tetra[®] System

an industry-renowned instrument for monitoring GPCRs and ion channels



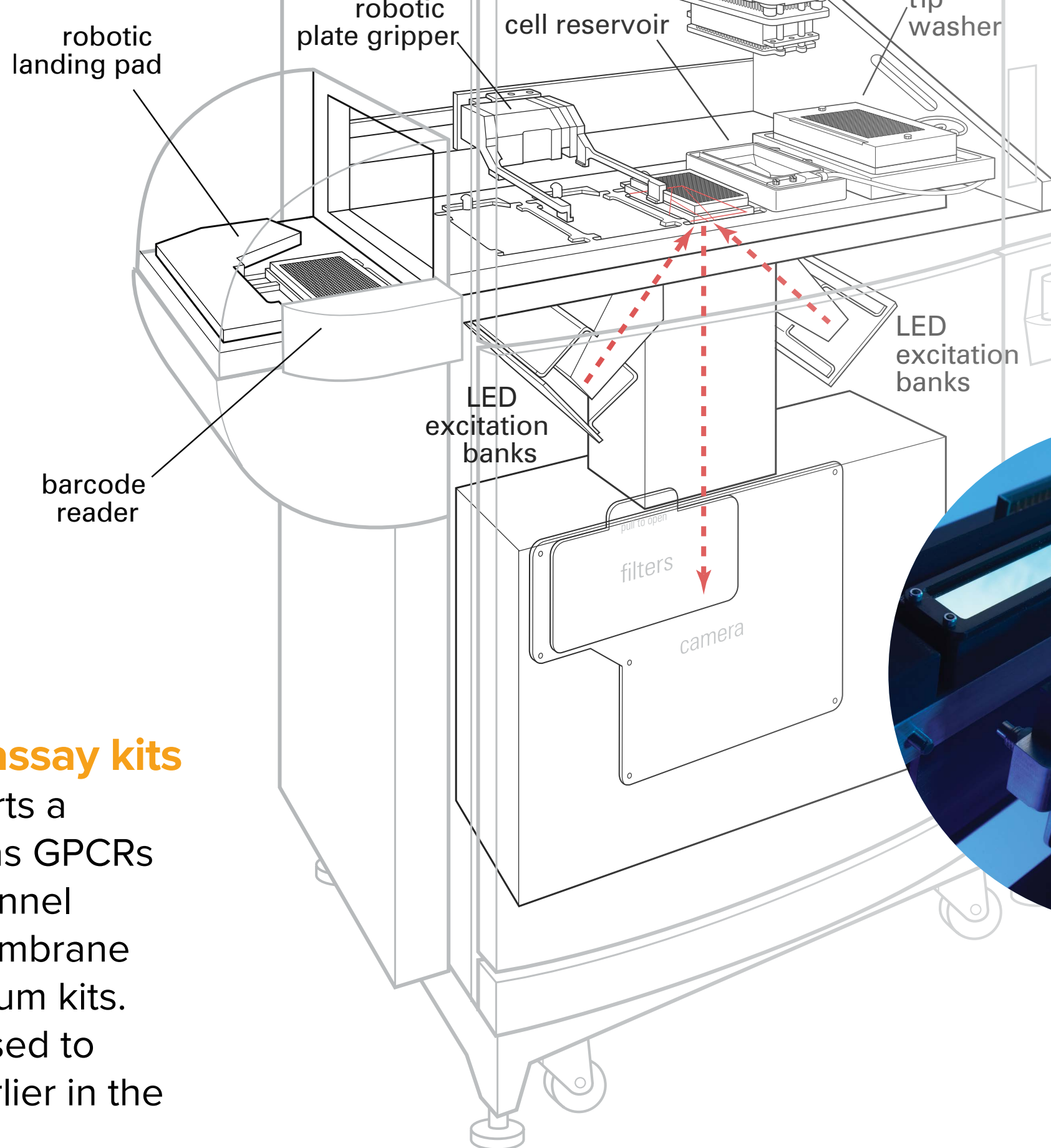
User-exchangeable 96-, 384-, and 1536-well pipettors

Install pipettors in minutes without tools. Pipetting accuracy and precision are assured with automatic pipettor identification and alignment.



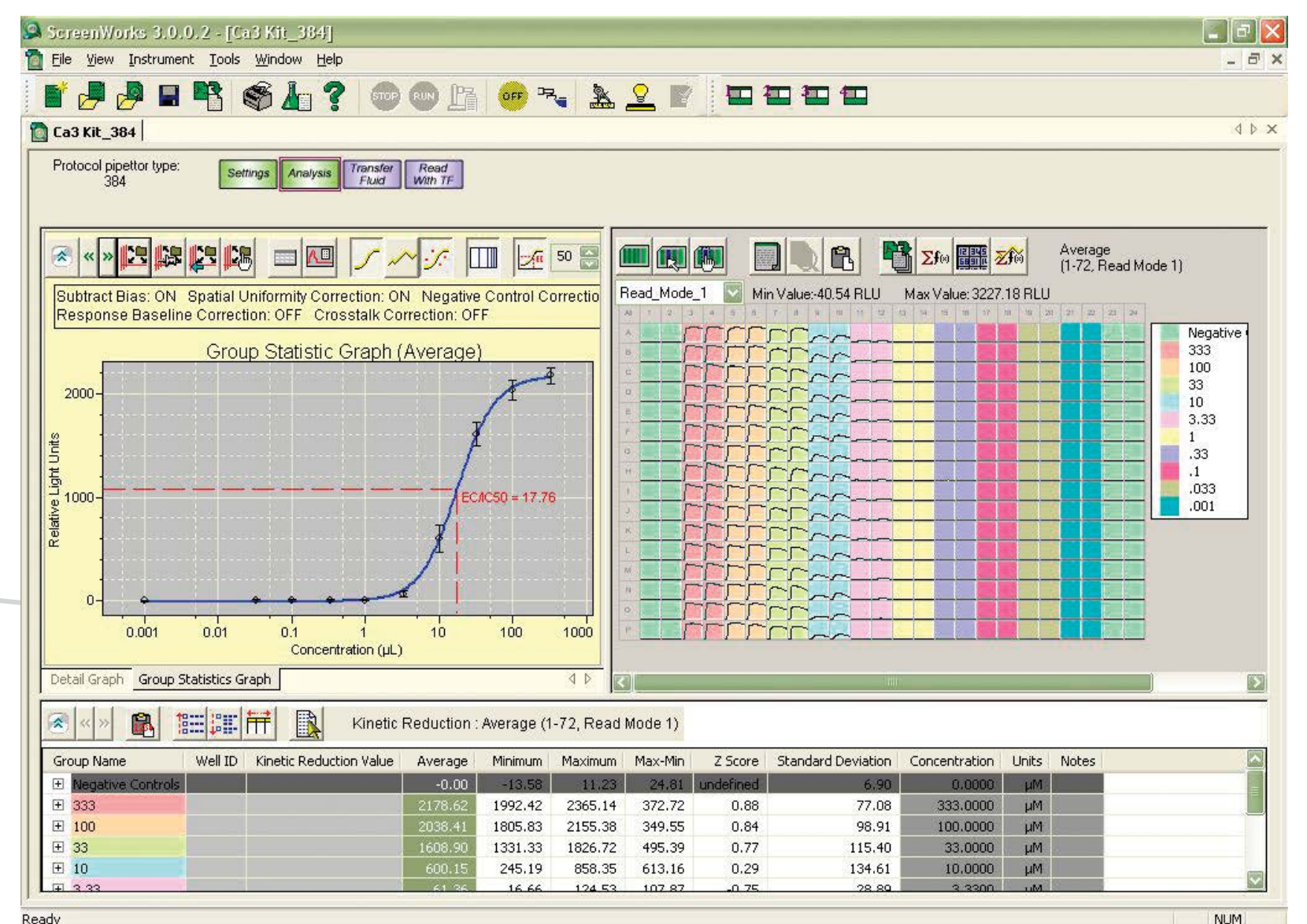
Fast, effective high-throughput cellular screening

Simultaneous liquid transfer to approximately 2,000 cells per well in a 1536-well plate is achieved using a proprietary, contact-based elastomeric technology. Further optimize speed with our Advanced Workflow Engineering Solutions team.



Supports a wide range of assay kits

The FLIPR Tetra System supports a wide range of assay kits such as GPCRs using Calcium kits and Ion Channel assays using optogenetics, Membrane Potential, Calcium, and Potassium kits. Cardiomyocytes can also be used to evaluate compound effects earlier in the drug discovery process.



Intuitive, user-friendly software interface

Easily create new protocols and access data. Analyze data in ScreenWorks software and create EC₅₀ or IC₅₀ curves with just a few clicks or set up auto export option for third party analysis.

Exclusive, configurable optics

Two camera options—standard for fluorescence only, or aequorin for both fluorescence and luminescence—allow you to tailor the system to your needs.

Click on the respective assay header for more information

GPCR	Membrane Potential	Cardiomyocytes	Potassium	Ligand-gated Calcium Channels	Optogenetics	Aequorin
Monitor live-cell calcium mobilization by GPCRs and ion channels	Observe real-time membrane potential changes associated with ion channel activation and ion transporter proteins	Observe calcium flux to assess cardiotoxicity in the early stages of drug discovery	Measure functional activity of ligand- and voltage-gated potassium channels	Measure the change of global intracellular calcium, in response to the activation of ion channels	Control membrane voltage in a reversible and precise fashion for screening state-dependent calcium channel blockers	Measure calcium flux using a luminescence assay for identifying early GPCR and ion channel leads during drug discovery

For more information, visit us at moleculardevices.com/FLIPR