

Ion Channel Retreat

Exploitation of Barracuda Automated Electrophysiology Platform to Improve Assessment of Cardiac Liability

Matt Bridgland-Taylor, Discovery Sciences

AstraZeneca R&D Alderley Park, UK

June 2013

matthew.bridgland-taylor@astrazeneca.com



Diversity of drug-induced cardiovascular toxicity

- Major cause of drug attrition in pre-clinical and clinical development
- Cause of serious adverse drug reactions (ADRs) and withdrawal from the market.

Heart

• Functional

- Arrhythmia
- Contractility

QT / QRS/ PR
prolongation

Cisapride

Blood Vessels

• Functional

- Blood pressure changes
- Orthostatic hypotension
- Oedema

• Structural

- Remodelling
- Vasculopathies

Cardiomyopathy
Myocardial infarction

Heart failure

Heart

• Structural

- Necrosis
- Fibrosis
- Hypertrophy
- Valvulopathy

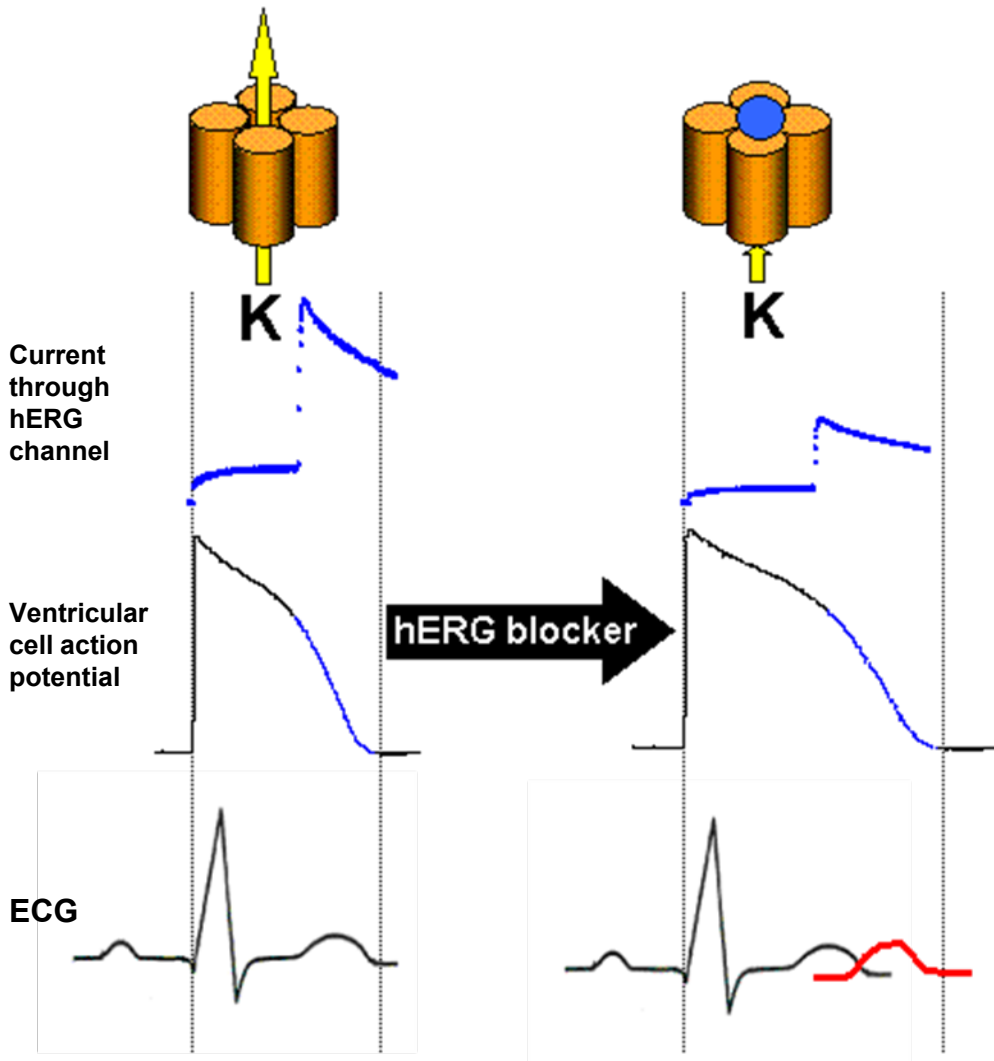
Blood

- Thrombosis
- Platelet function
- Anaemia/ RBC haemolysis
- Immune system



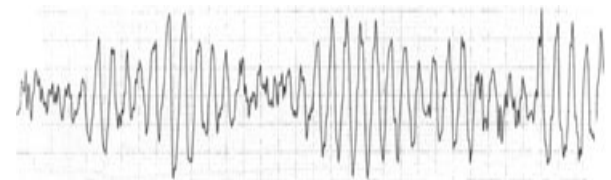
Drug-induced QT interval prolongation

Biological background



- hERG channel block prolongs ventricular cell action potentials
- Seen as QT interval prolongation on ECG
- QT interval prolongation is linked to proarrhythmic risk (Torsades de Pointes)
- hERG block is most common mechanism
- QT interval has to be corrected for heart rate changes (QTc)
- If drug significantly increases heart rate, this can lead to an apparent QTc increase

Torsades de Pointes (TdP)

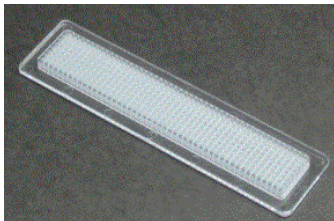
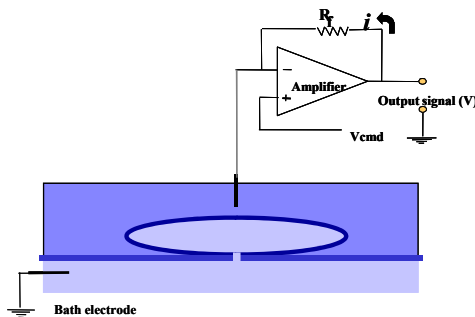


IonWorks versus GLP hERG

Technical differences

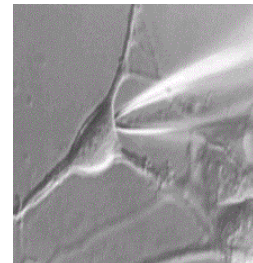
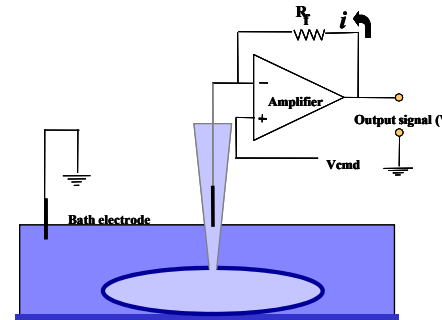
- IonWorks

- 384-well, plate-based
- Plastic surface area:buffer volume ratio, high
- Intermittent voltage clamp
- Rate of onset of block cannot be measured*
- No measurement of drug concentration
- Therefore, potency based on nominal concentrations



- Good Lab Practice (GLP)

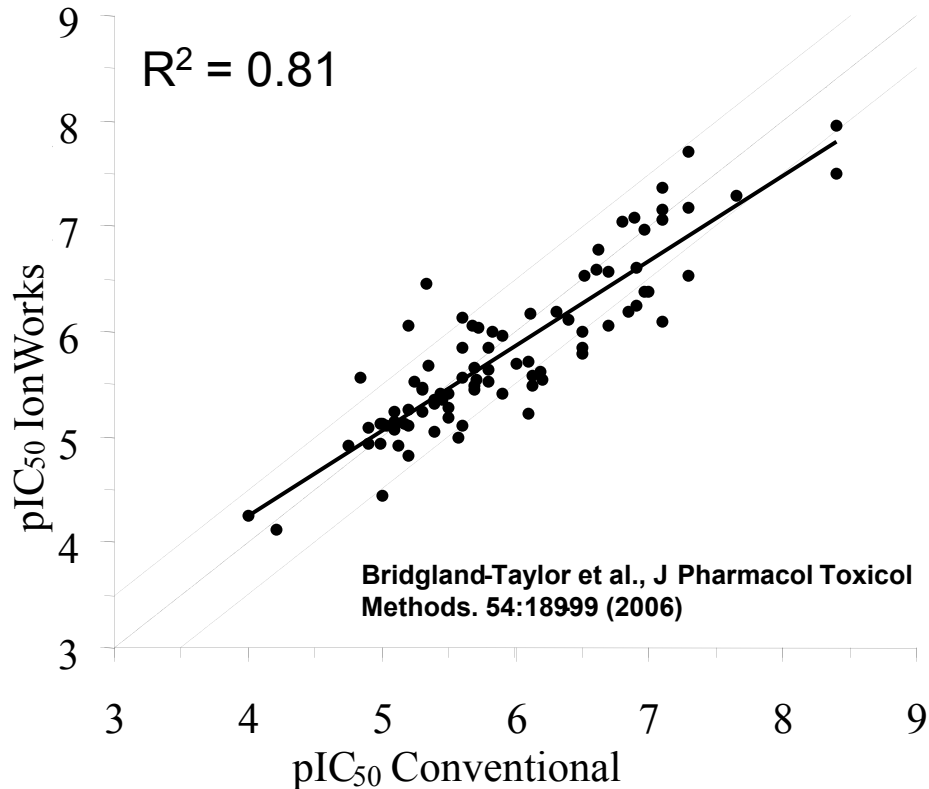
- Microelectrode-based conventional patch clamp
- Plastic surface area:buffer volume ratio, low
- Continual voltage clamp
- Rate of onset of block can be measured
- Drug concentration measured
- Therefore potency based on measured concentrations



GLP versus IonWorks hERG

Data correlations

- Initial IonWorks validation prior to screen launch in 2003
 - For 89 compounds comparing IonWorks to microelectrode method, correlation was generally good with a few exceptions (for example AZ1)



Potential explanations for a difference in potency between IonWorks and GLP hERG are:

- The actual concentration in IonWorks is less than the nominal value
- The compound is a very slow on-set blocker such that its effect is underestimated in IonWorks
- The lack of continuous voltage clamp in IonWorks

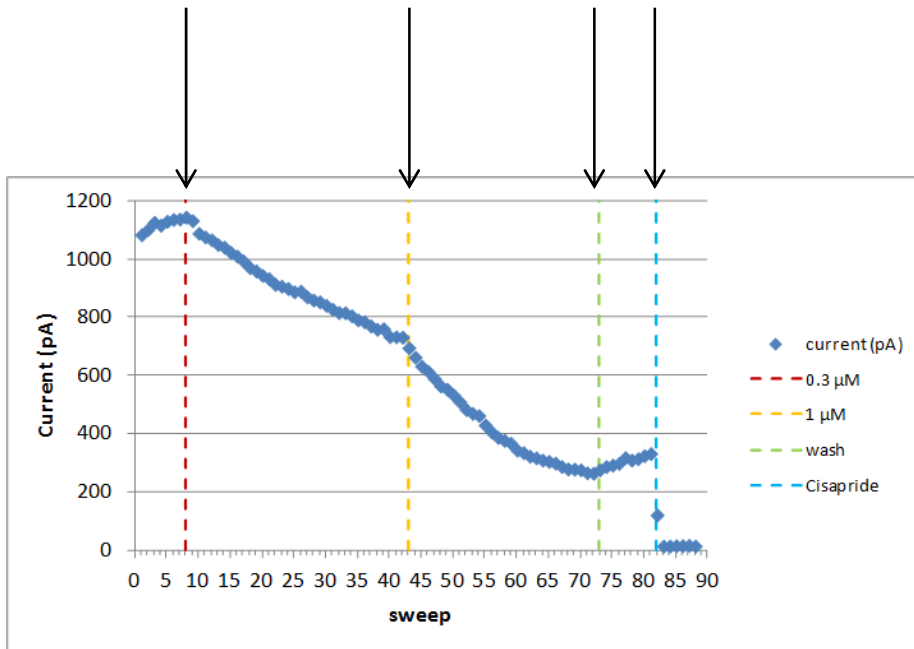


AZ1 IonWorks:GLP hERG data

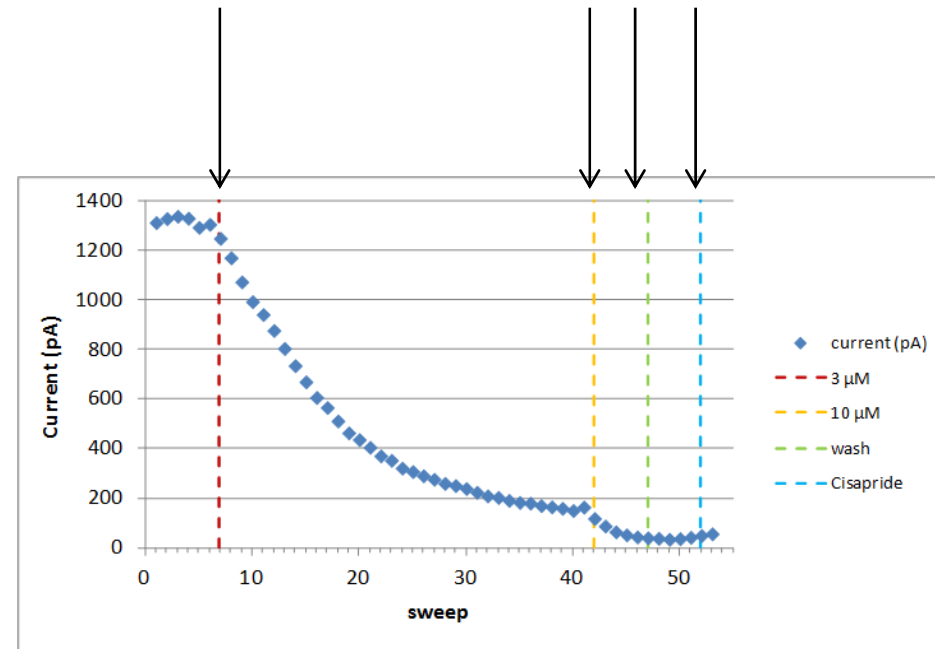
IC₅₀ of 22.1 μM (n=6) for standard IonWorks-based assay (200 sec incubation)
Compared to IC₅₀ of 0.25 μM in GLP-hERG (n=4) – is it a slow onset compound?

Trace profile characteristic of slow onset compound

Compound additions



Compound additions

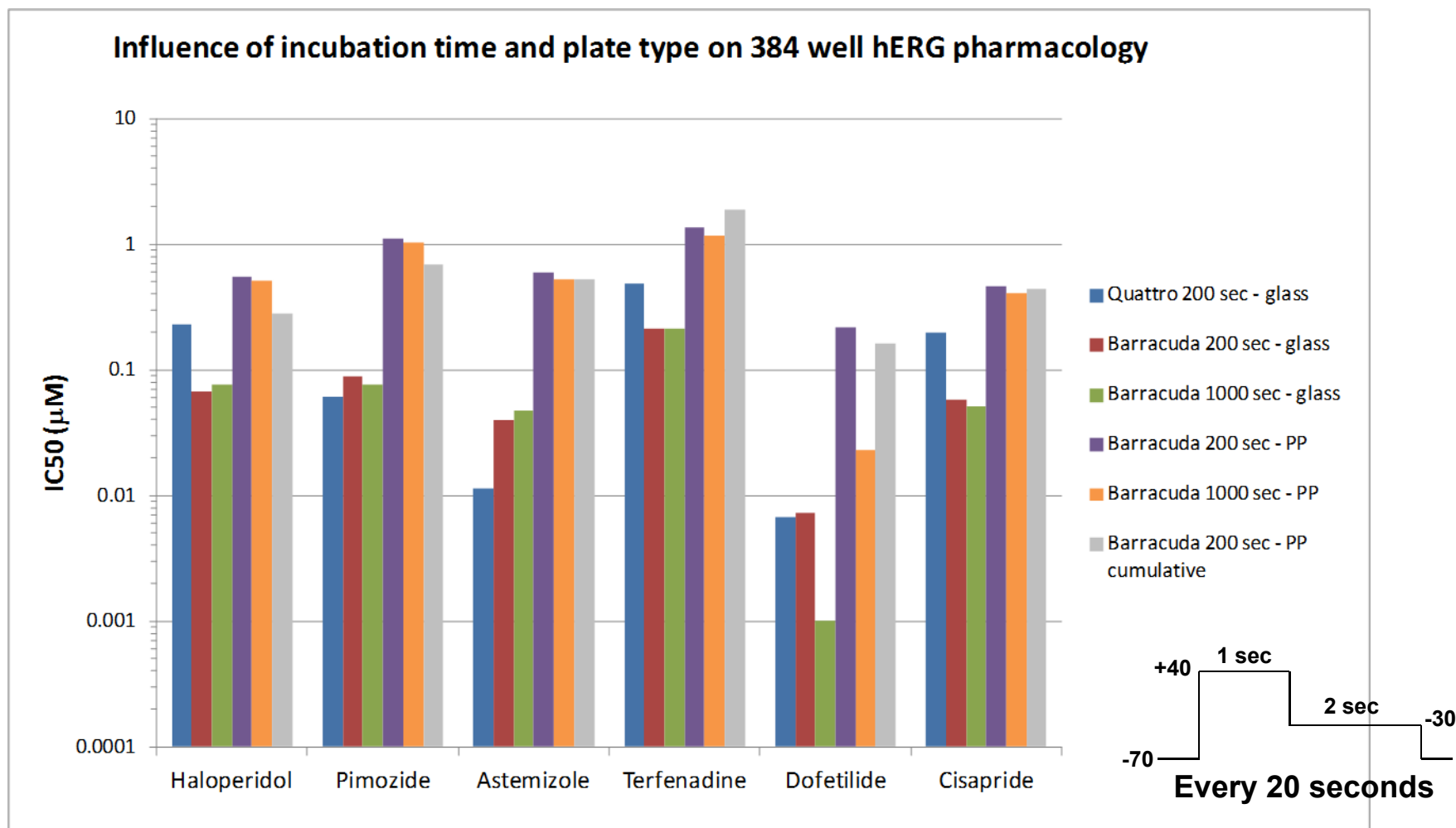


The next generation IonWorks device (Barracuda), like the GLP hERG assay, enables continuous voltage clamp and measurement of on-set of block



Barracuda validation - comparing incubation time

Assessing compound plate type and incubation time of hERG standards



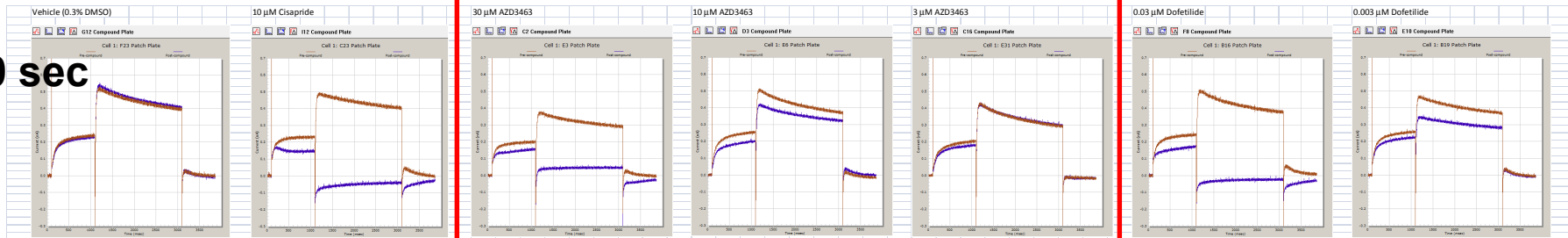
Is AZ1 similar to dofetilide (slow onset blocker)?



Side by side IonWorks vs Barracuda traces

IonWorks

200 sec

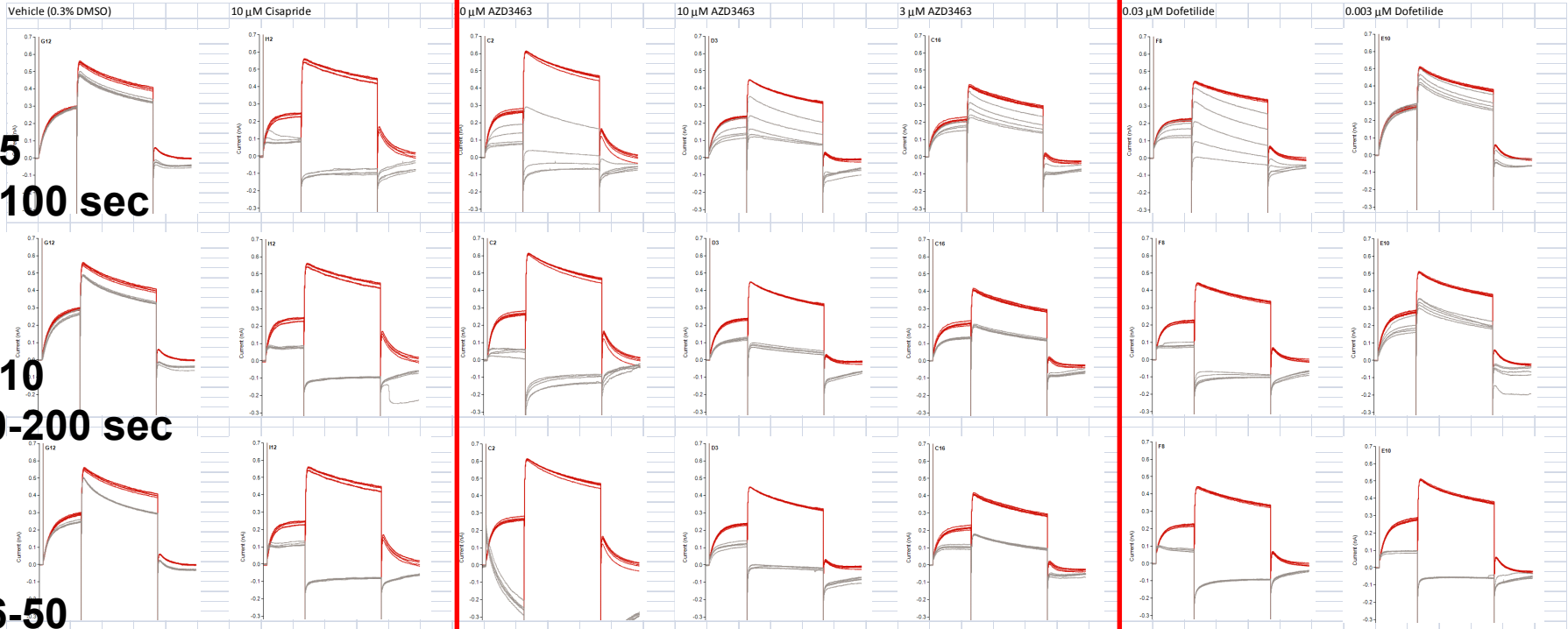


Barracuda

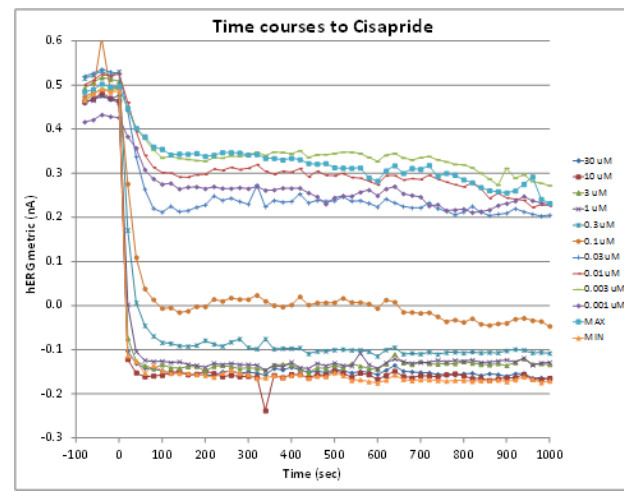
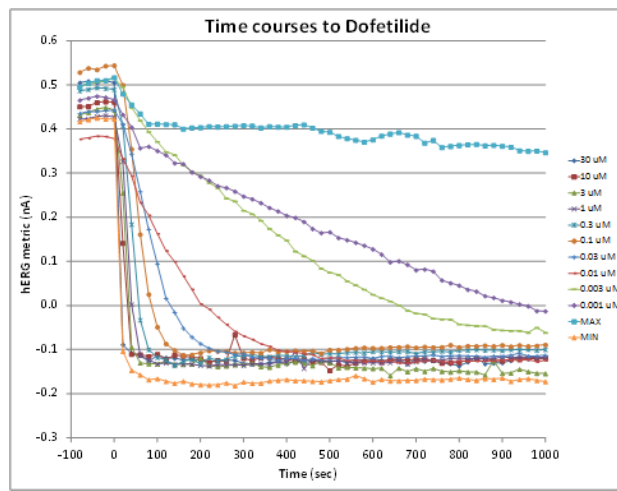
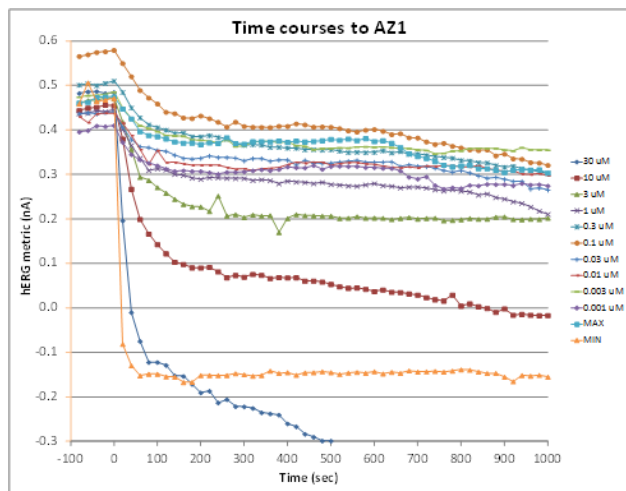
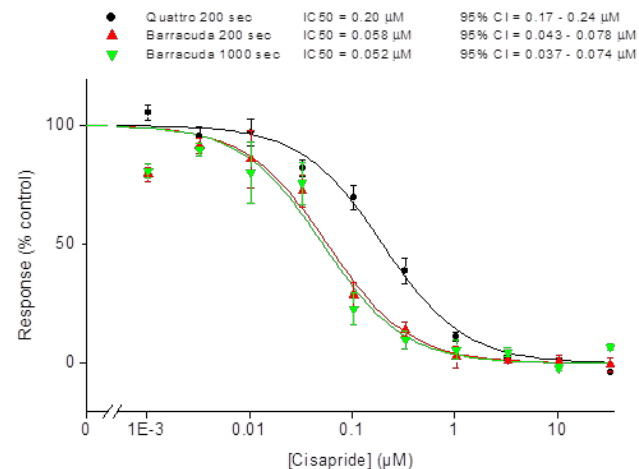
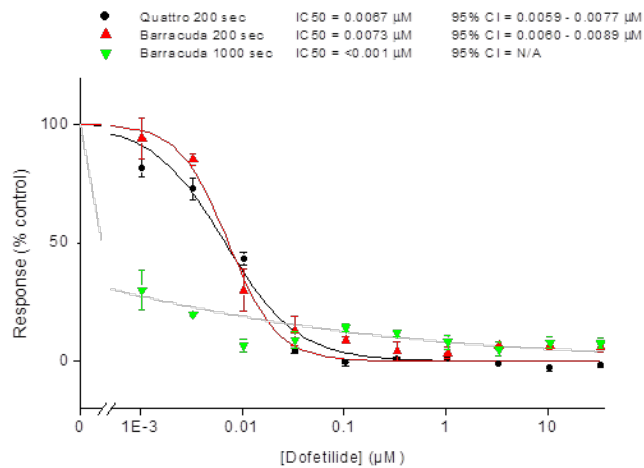
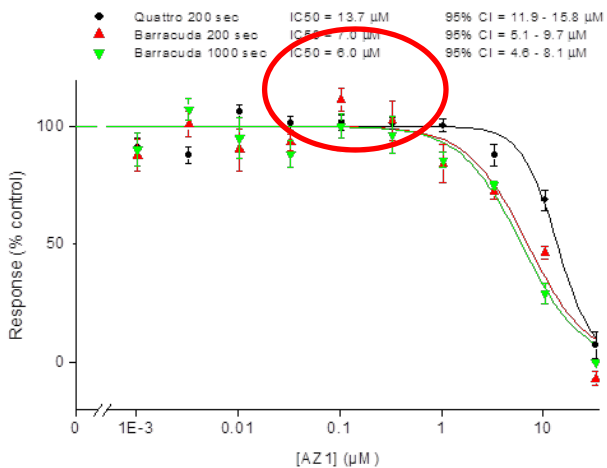
s1-5
20-100 sec

s6-10
120-200 sec

s46-50
920-100 sec



Data using glass-coated compound plates (n=4)



Comparing GLP, IonWorks & Barracuda

Summary of AZ1 hERG IC₅₀ values using different methods

	IonWorks in HT mode: 96 well polypropylene (200 sec)	IonWorks in PPC mode: 384 well glass coated compound plate (200 sec)	GLP-hERG: minimal amount of plastic	Barracuda in PPC mode: 384 well glass coated compound plate (1000 sec)
Based on nominal concentration	22.1 µM	13.7 µM	0.42 µM	6.0 µM (7.0 µM at 200 sec plus detection of slow on-set of block)
Based on measured concentration	5.3 µM (extraction of contents from PatchPlate for DMPK analysis)	Not Determined	0.25 µM	**Not Determined

**** = if assume similar cell concentration & plastic surface area between IonWorks & Barracuda, then IC₅₀ should drop to ~1.4 µM (5.6 fold from GLP-measured).**



Summary of work with AZ1

- **IonWorks Barracuda™ was able to detect the very slow on-set of hERG block when testing AZ1. Some of the remaining difference in potency when compared to the GLP-hERG data is due to the compound sticking to the PatchPlate substrate or the difference in cell concentration.**
- **If we assume a similar drop in potency seen on IWQ when using the IWB PatchPlate, the potency of AZ1 is only 5.6 fold greater than the GLP-hERG IC₅₀ value (~1.4 versus 0.25 μM).**
- **Gillie et al (2013) have reported the effects of longer exposure to Dofetilide in the IonWorks Barracuda™ system and here we have also demonstrated it is a very slow on-set blocker of the hERG channel. We have also seen that Dofetilide potency is influenced by increasing the incubation time with the IonWorks Quattro system (data not shown).**
- **Using IonWorks Barracuda™ for routine hERG screening is delivering a better assessment of compound block at this important ion channel.**



Confidentiality Notice

This file is private and may contain confidential and proprietary information. If you have received this file in error, please notify us and remove it from your system and note that you must not copy, distribute or take any action in reliance on it. Any unauthorized use or disclosure of the contents of this file is not permitted and may be unlawful. AstraZeneca PLC, 2 Kingdom Street, London, W2 6BD, UK, T: +44(0)20 7604 8000, F: +44 (0)20 7604 8151, www.astrazeneca.com

