

Introduction

Drug development costs exceed \$172m and increase to over half a billion dollars when clinical trial failures are taken into account¹. It can take up to 20 years to successfully bring a drug to market, so it is critical to identify any safety liabilities early in the drug discovery process to ensure valuable time and money are not invested unnecessarily in time-consuming and expensive preclinical animal and clinical trials.

Despite the importance of identifying safety liabilities early in drug discovery, over the past 2 decades, the proportion of venture capital investments in early stage (including in vitro safety panels) has decreased from **50%** (2006) to **16%** (2023) while increasing for later stage (including clinical trials) from **20%** to **54%**^{2,3}. This indicates an understandable trend towards expensive, **high-risk** but potentially lucrative, clinical trials.

The goal of this study is to demonstrate that early-stage in vitro safety panels are still critical steps in the drug development process both fiscally and scientifically as these identify off-target effects of drug candidates before advancing to expensive and time-consuming animal and human testing.

Methods

- 20 drug candidates from failed oncology clinical trials (2001-2025)⁶
 - Reason for termination: **severe adverse effects**
- Compounds screened at 10 μM against 44 safety liability targets in InVEST44™ and 70 kinases in Diversify™
- Table 1 lists the targets for **InVEST44™ in green**.

Target families include:

- G-protein coupled receptors (GPCRs)
- Ion channels
- Transporters
- Nuclear receptors
- Kinases
- Other enzymes

Assay types include:

- Competitive radioligand binding
- Fluorescence polarization
- Manual patch electrophysiology
- FLIPR calcium mobilization
- Cell reporter
- Enzymatic activity
- ADP-Glo

InVEST™ In Vitro Safety Suite

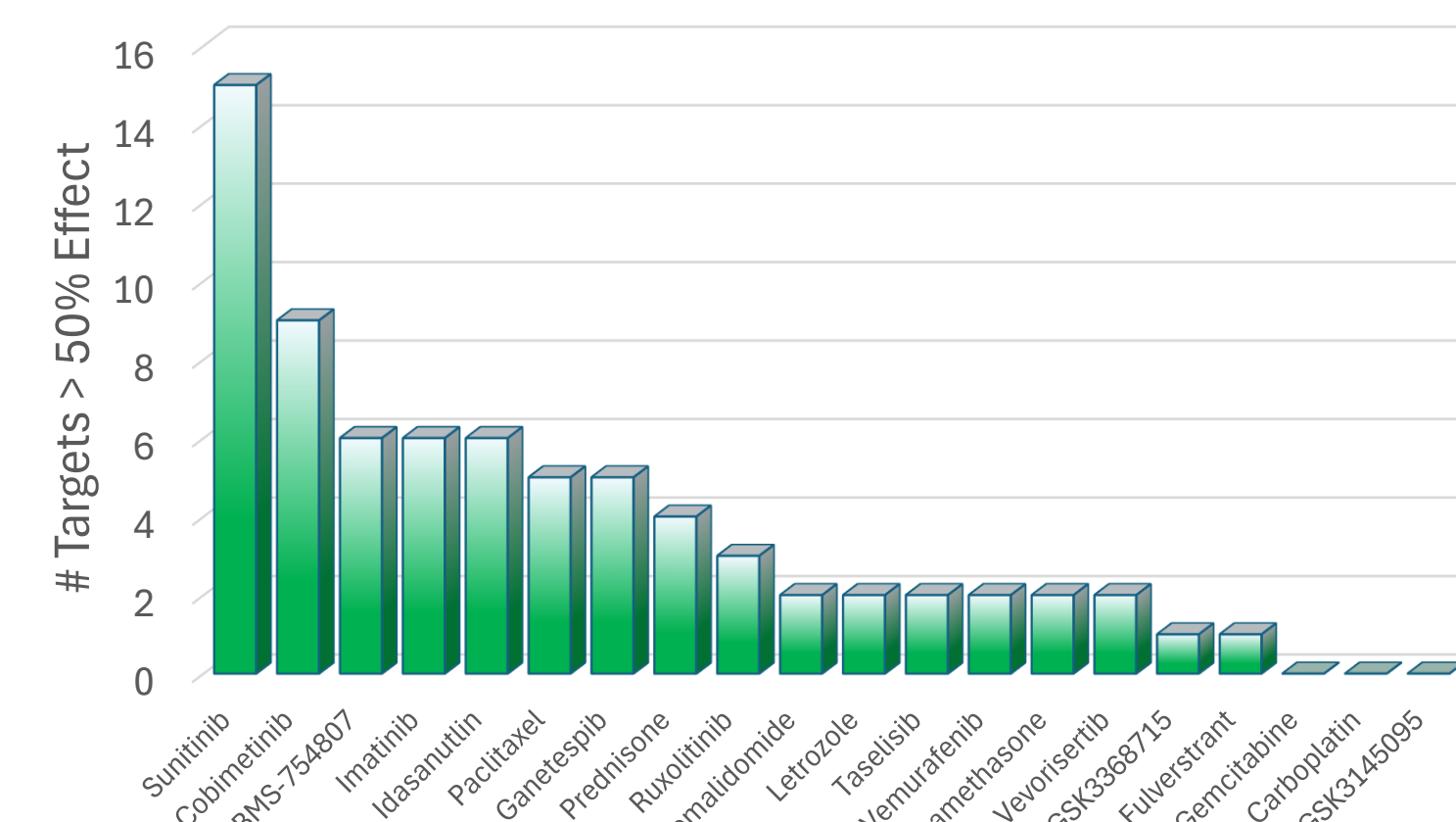
At Reaction Biology we offer a suite of in vitro safety panels – **InVEST18™**, **InVEST44™**, **InVEST59™** and **InVEST77™** - providing an opportunity to identify potential toxicities of lead compounds in a cost-effective manner and at an early stage in drug development.

The targets selected for **InVEST18™**, **InVEST44™**, **InVEST59™** and **InVEST77™** are based on surveys across 4 major pharmaceutical companies⁴ and 18 companies overseen by IQ DruSafe⁵. These targets are crucial in both core systems (CNS, CV, respiratory) and secondary organs (GI, renal).

Results

Failed clinical trial candidates exhibit off-target safety liabilities in InVEST44™ panel

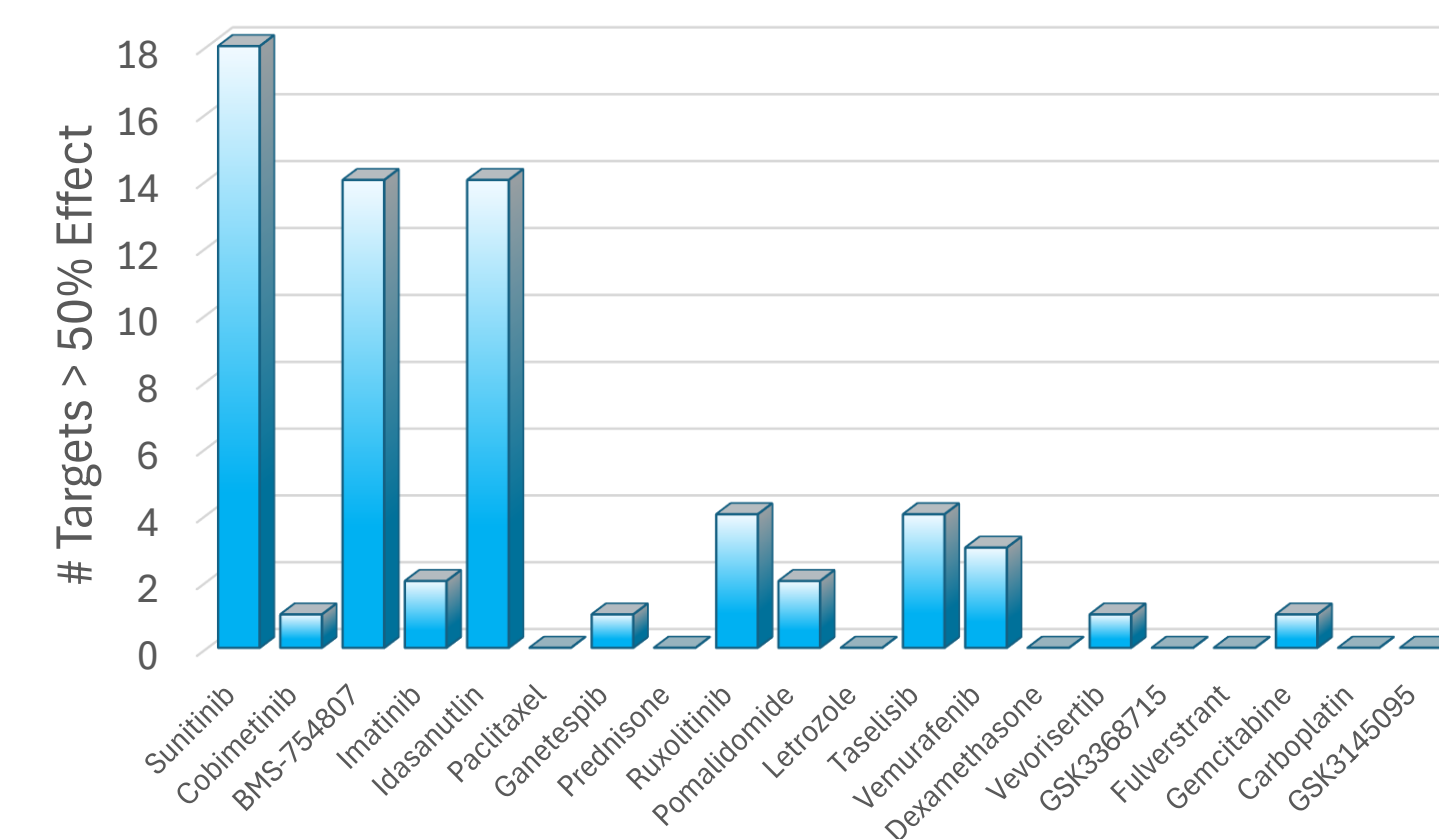
Off-Target Effects Identified by InVEST44™



17 of the 20 drugs elicited >50% effect on at least 1 target in the InVEST44™ panel, including 8 with 4+ target interactions. One drug candidate interacted significantly with 15 safety targets.

Failed clinical trial candidates exhibit multiple kinase interaction in kinase Diversify™ panel

Effects of Kinases Identified Diversify™



12 of the 20 drugs elicited >50% effect on at least 1 target in the Diversify™ panel, including 5 with 4+ target interactions. One drug candidate interacted significantly with 18 kinase targets. The Diversify™ panel comprises 70 kinases selected to represent multiple kinase families. For a complete list of targets, see <https://www.reactionbiology.com/list-kinase-targets-us-facility/?panel=Diversify%20Panel>

InVEST™ Targets

Target Family	InVEST18	InVEST44	InVEST59	InVEST77	
Kinase			ATM	ATM	
			AURKA	AURKA	
			AURKB	AURKB	
			BRD4	BRD4	
			EGFR	EGFR	
			FGFR1	FGFR1	
			FLT1	FLT1	
			GSK3B	GSK3B	
			INSR	INSR	
			KDR	KDR	
			KIT	KIT	
		Lck	Lck	Lck	
			MAP2K7	MAP2K7	
			PDK1	PDK1	
			PIK3CG	PIK3CG	
			PTK2	PTK2	
			ROCK1	ROCK1	
			ROCK2	ROCK2	
		SRC	SRC		
		STK35	STK35		
GPCR		ADORA2A	ADORA2A	ADORA2A	
		ADRA1A	ADRA1A	ADRA1A	
		ADRA2A	ADRA2A	ADRA2A	
		ADRB1	ADRB1	ADRB1	
		ADRB2	ADRB2	ADRB2	
		CNR1	CNR1	CNR1	
			CNR2	CNR2	
			CCK1	CCK1	
			CCK2	CCK2	
		D1	D1	D1	
			D2	D2	
			D3	D3	
			D3	D3	
			ETA	ETA	
			GPBAR1	GPBAR1	
		H1	H1	H1	
			H2	H2	
			OPRD1	OPRD1	
		OPRK1	OPRK1		
		OPRM1	OPRM1		
		OPRM1	OPRM1		
		CHRM1	CHRM1		
		CHRM2	CHRM2		
Ion Channel					
		CaV1.2	CaV1.2	CaV1.2	
		GABR	GABR	GABR	
			HTR3	HTR3	
			HTR3	HTR3	
		KV11.1	KV11.1	KV11.1	
			CHRNA1	CHRNA1	
			CHRNA4	CHRNA4	
			NMDA	NMDA	
		SCN5A	SCN5A	SCN5A	
	Enzyme				
			ACHE	ACHE	ACHE
				COX1	COX1
				COX2	COX2
				CTSD	CTSD
				MALT1	MALT1
				MAOA	MAOA
				MAOB	MAOB
			PDE3A	PDE3A	
			PDE4D2	PDE4D2	
			PDE4D2	PDE4D2	
		DAT	DAT	DAT	
			NET	NET	
			SERT	SERT	
			AR	AR	
			ESR1	ESR1	
			GR	GR	
			PGR	PGR	
		RARA	RARA		
Nuclear Receptor					

Table 1: Safety liability targets for InVEST18™, InVEST44™, InVEST59™ and InVEST77™

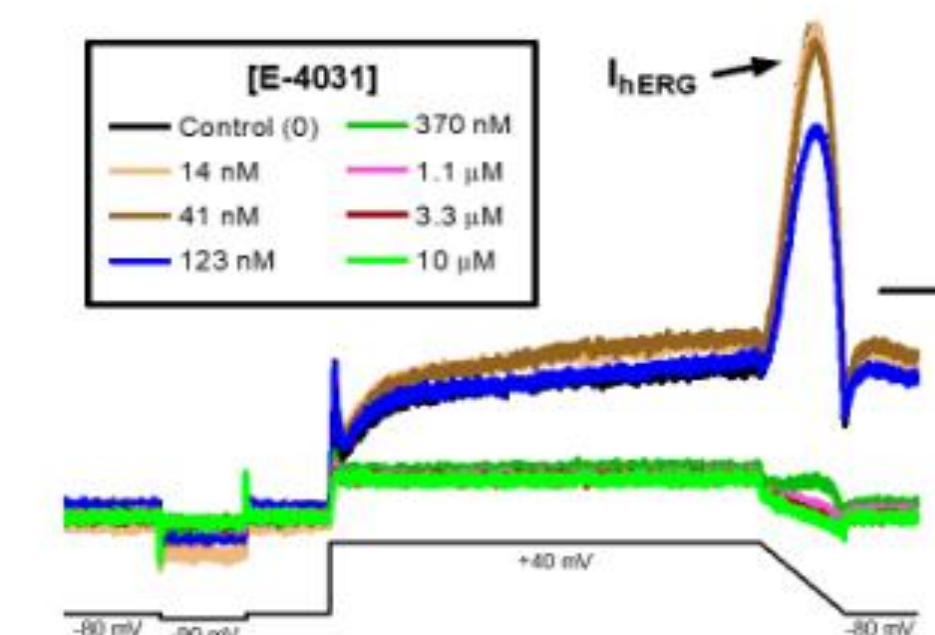
Conclusion

In vitro safety panels, such as **InVEST44™**, are critical in identifying off-target effects of lead compounds prior to time-consuming, expensive animal testing and failed clinical trials.

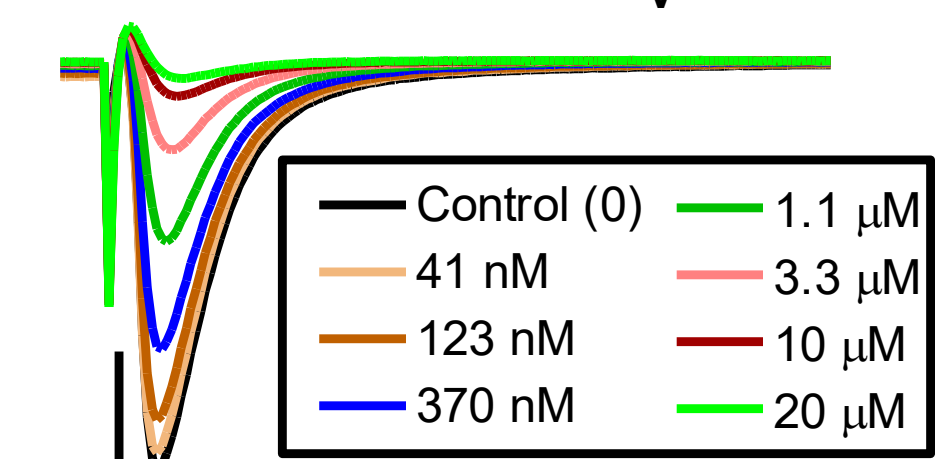
By providing cost-efficient, early-stage identification of off-target effects by lead compounds, in vitro safety panels, such as **InVEST44™**, have become integral components of drug discovery and warrant increased early-stage consideration and investment in the future.

Other In Vitro Safety Offerings

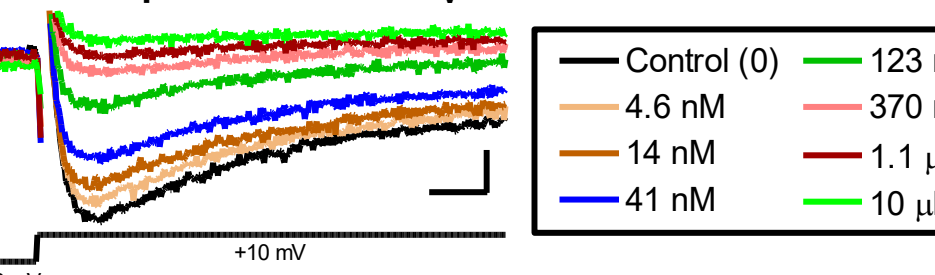
Reaction Biology offers manual and automated patch whole-cell electrophysiology for **cardiac safety**, including **hERG**, **NaV1.5** and **CaV1.2**. Ask for more details.



TTX Block of NaV1.5



Nifedipine Block of CaV1.2



Scale bars: hERG: 100pA, 1s; NaV1.5: 5nA, 1ms; CaV1.2: 100pA, 5ms

References

- Sertkaya et al. (2024) JAMA Netw Open 7(6):e415445. DOI: 10.1001/jamanetworkopen.2024.15445.
- BIO. (2015). *The Biopharma Industry Investment Report 2015*. Biotechnology Innovation Organization
- Larka, M. (2024). *Trends in Venture Capital Funding for Biopharma Startups*. GlobalData
- Bowes et al. (2012) Nat Rev Drug Discov 11(12):909-22. DOI: 10.1038/nrd3845.
- Brennan et al. (2024) Nat Rev Drug Discov 23(7):525-545. DOI: 10.1038/s41573-024-00942-3.
- clinicaltrials.org

Contact Information

Julian Woollorton, PhD
Director, In Vitro Safety

Reaction Biology Malvern
1 Great Valley Parkway,
Suite 2
Malvern, PA 19355

+1 610-722-0247
Julian.Woollorton@reactionbiology.com
www.reactionbiology.com