

MAKE TIME FOR THINGS
THAT MATTER

ELISA automation
with Crocodile
saves you time
and money
















CALCULATE YOUR ELISA ROI

Learn how you can save precious time and money even when performing low throughput ELISA

ELISA protocols are tedious and involve several time-consuming manual steps. These take up valuable time of the staff, that could be used more productively.

The Crocodile miniWorkstation provides a simple and affordable automation solution for low throughput ELISA, freeing up valuable human resources for other important tasks.

TYPICAL PROTOCOL	MANUAL OPERATION	WITH CROCODILE	Cost factors	
 Setup			How many plates (whole plate or part of a plate) do you process per day (total plates of typical ELISA or similar assays)?	<input type="text"/>
 Prepare	Operator busy	Operator busy	How much is the hourly rate of the scientist or technician typically running your assays (full costs)?	<input type="text"/>
	Walk/Wait	Walk/Wait		
 Incubate	Operator free			
	Walk/Wait			
 Wash	Operator busy			
 Dispense	Walk/Wait			
 Incubate	Operator free	Operator free		
	Walk/Wait			
 Wash	Operator busy			
 Dispense	Walk/Wait			
 Incubate	Operator free			
	Walk/Wait			
 Dispense				
 Measure	Operator busy	Walk/Wait		
 Tidy up		Operator busy		
 Shutdown				
			Washing	
			How much time do you spend in washing your plate in your typical assay run? (Total in minutes)	<input type="text"/>
			Dispensing	
			How much time do you spend in dispensing reagents in your typical assay run? (Total in minutes)	<input type="text"/>
			Walk / Wait / Transfer	
			How much time do you spend overall waiting that processing steps finish before you can continue with the next step in one typical assay run, not counting incubation times (for example, if time between steps is too short to perform any other tasks and you just wait)? (Total in minutes)	<input type="text"/>
			How much time do you spend overall walking between steps in one typical assay run (transferring your plate between instruments, walking between workplaces...)? (Total in minutes)	<input type="text"/>
			MANUAL OPERATION	WITH CROCODILE
OPERATOR TIME PER ASSAY [min]			<input type="text"/>	<input type="text"/>
OPERATOR TIME PER YEAR [h]			<input type="text"/>	<input type="text"/>
OPERATOR COSTS PER YEAR			<input type="text"/>	<input type="text"/>
TIME SAVINGS PER YEAR USING THE CROCODILE [h]				<input type="text"/>
COST SAVINGS PER YEAR USING THE CROCODILE				<input type="text"/>

Get a quote now

Contact our automation specialists

*We use the term “operator” for short, meaning the person performing the assay (lab technician, scientist...).

Steps of the protocol not entered in the calculator are assumed to need roughly the same time with or without the Crocodile (plate setup, incubation and measurement times...).

Operator costs and cost savings are calculated using the same currency under “Cost factors”.