

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

Setup

Prepare

Incubate

Wash

Dispense

Incubate

Wash

Dispense

Incubate

Dispense

Measure

Tidy up

Shutdown

MANUAL
OPERATION

Operator busy

Walk/Wait

Operator free

Walk/Wait

Operator busy

Walk/Wait

Operator free

Walk/Wait

Operator free

Walk/Wait

Operator busy

WITH
CROCODILE

Operator busy

Walk/Wait

Operator free

Walk/Wait

Operator busy

Cost factors

How many plates (whole plate or part of a plate) do you process per day (total plates of typical ELISA or similar assays)?

How much is the hourly rate of the scientist or technician typically running your assays (full costs)?

Washing

How much time do you spend in washing your plate in your typical assay run? (Total in minutes)

Dispensing

How much time do you spend in dispensing reagents in your typical assay run? (Total in minutes)

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)

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)

MANUAL
OPERATION

WITH
CROCODILE

OPERATOR TIME PER ASSAY [min]

OPERATOR TIME PER YEAR [h]

OPERATOR COSTS PER YEAR

TIME SAVINGS PER YEAR USING THE CROCODILE [h]

COST SAVINGS PER YEAR USING THE CROCODILE

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”.

For Research Use Only. Not for use in diagnostic procedures

© 2023 Berthold Technologies. All rights reserved. All trademarks are the property of Berthold Technologies and its subsidiaries unless otherwise specified.