

Example data for the Crocodile ROI Calculator

To simplify filling in the data in the Crocodile ROI Calculator, we provide 3 pre-filled examples with the timing data as estimated by 3 different customers.

IMPORTANT: hourly rates might be very different depending on country, position of the person performing the ELISA, type of company/institution, and so on. Please try to input a rate that is representative for you (considering how much the hour costs to the company/institution, not how much the person earns).

The 3 cases are:

1. In case 1, both washing and reagent dispensing are performed using a multipipette or similar repeating device. During incubations, the operator works close to where the ELISA is being performed, so not much time is wasted walking between workplaces. However, combining many different activities during incubations makes planning difficult, so some time is wasted because sometimes the time available between activities is too short to do anything meaningful.
2. In case 2 reagent dispensing is performed with a single-channel pipette, and washing with a multi-channel pipette. During incubations, the person works mainly in the office, and can plan other activities with a minimal waste of time. However, the office is in a different floor, so they spend a significant amount of time walking from the lab to the office and back.
3. In case 3, reagents are dispensed using a multipipette or similar repeating device, and washing is performed using a wash bottle, resulting in reduced hands-on time. However, the last incubation is typically 10 minutes only, which makes it very difficult to use it to perform other meaningful tasks. Thus, this time is added to the time spent waiting.

If you want a quick estimation, just pick the case that is more similar to your situation and enter your typical number of plates per day and hourly rate.

If you want to try to fill everything by yourself, but are unsure about the time you need for each step, use the data provided by our customers as reference:

Dispensing of reagents:

1. If using a multipipette or similar repeating device, 4-8 minutes.
2. If using a single channel pipette, 20-30 minutes.

Washing:

1. If using a multichannel pipette or wash bottle, 5-10 minutes.
2. If using a multipipette or similar repeating device, around 15 minutes.

Time waiting:

1. If using incubation times mainly working on the computer, 1-5 minutes.
2. If combining different experimental tasks, 5-10 minutes.
3. If the incubation with the substrate in your typical ELISA is short (10-15 minutes), it's possible that you can't use this incubation efficiently to perform other tasks. If this is your case, consider adding it to the time spent waiting.

Time walking:

1. If working during incubations in a workplace close to where the ELISA is running, 1-3 minutes.
2. If working during incubations in a different floor or building, 5-10 minutes or more (depending on the walking distance).