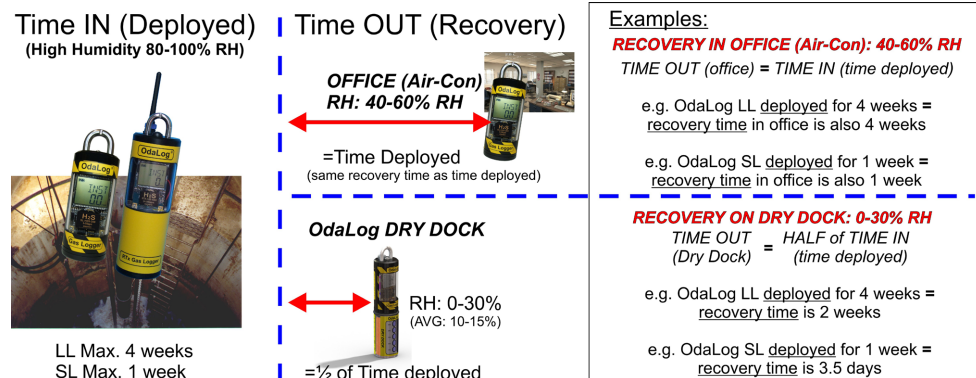


'TIME IN' - 'TIME OUT' DEPLOYMENT / RECOVERY

Due to the deployment / recovery time being inversely relational (each sensor recovers at the same 'rate' it absorbs moisture), the calculation for how long to 'dry' excess moisture from the sensor remains the same between both SL and LL model OdaLog Instruments. Use of the OdaLog Dry Dock, reduces the recovery time by half:



RECOVERING SATURATED DESSICANT BEADS

When the desiccant beads reach their limit of absorption, the colour will change to a deep green/blue. Often only the top layer of beads will change colour - in this case, replacing the rubber cover and allowing the beads to stand for 1-2 days equalise the moisture within the dry dock, making it ready for further use. Reactivation should be very infrequent (only small amounts of moisture in the sensor).

If reactivation is required, follow these steps: (1) Remove top cap from the Dry Dock; (2) Pour beads onto an oven tray, spreading evenly over tray surface; (3) Place tray into a fan-forced oven at 105-110°C until beads return to orange in colour; (4) Remove tray from oven and allow to cool for at least 5-10 minutes; (5) when safe to handle, pour beads back into Dry Dock and replace top cap and rubber cap. (**Note: beads will be extremely hot, and steam may be present - take all precautions.**)

FOR MORE INFO:

Contact your Authorised Distributor for more details
or email: sales@app-tek.com

ODALOG DRY DOCK QUICK USER GUIDE

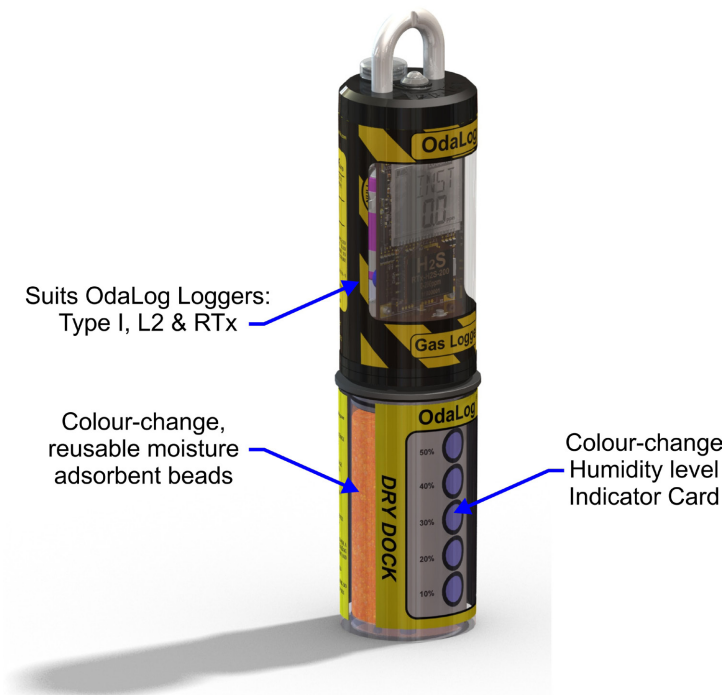
Package Contents:

OdaLog® Dry Dock (PN: 25-0391), Consisting of:

- OdaLog Dry Dock - Suits OdaLog Type I, L2 and RTx
- Highly-absorbent reusable colour-change desiccant beads (Orange-Blue, 200g)
- Humidity indicator card (indicates humidity levels between 10%-50%)
- Moulded (clear) Cover (keep fitted when Dry dock not in use)
- Quick User Guide (this document)

Accessories: SOLD SEPERATELY

- 12-0066, Replacement colour-change desiccant beads (Orange-Blue, 200g)



WARNINGS

- **'Over-drying' (leaving instrument on Dry Dock too long) may cause damage to the sensor** (affects performance), do not dry for longer periods than recommended.
- Do not attempt to 'recover' (heat in oven / microwave) desiccant beads while in the OdaLog Dry Dock - **beads must be removed from the plastic body before recovering / heating**: plastic parts are NOT microwave-safe.

GENERAL DISCLAIMER

The *OdaLog® Dry Dock* is sold subject to our standard terms and conditions of sale and any written warranty given by App-Tek. App-Tek does not accept any other liability either directly or indirectly for any losses suffered in connection with the use and application of the *OdaLog® Dry Dock*, whether or not in accordance with any advice, specification, recommendation or information given by it or as a result of any errors or omissions in this *Quick User Guide*.

USING THE ODALOG DRY DOCK

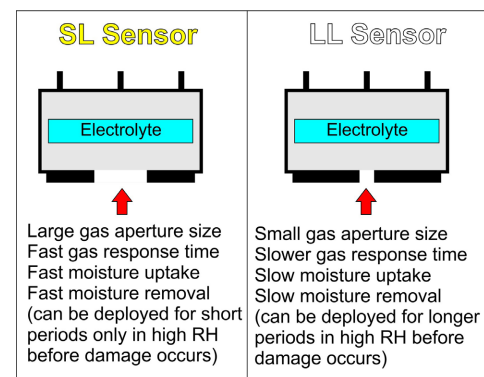
1. Check the humidity level inside the OdaLog Dry Dock is low: beads should be orange in colour, and at least the 50/40/30 RH circles on the indicator card must be blue.
2. Check the OdaLog sensor filter/grill for obstructions (replace if clogged).
3. Remove the rubber cover from the dry dock, and gently place OdaLog onto the Dry Dock.
4. Calculate the time required for recovery, and ensure that the OdaLog does not exceed this period (over-drying will affect sensor performance). Use a countdown timer if available. (See 'Time in' - 'Time Out' Deployment / Recovery' section on following page)
5. Once the recovery time is complete, remove the OdaLog from the Dry Dock. Immediately replace the rubber cap to the Dry Dock (keeps beads from absorbing humidity from air).

Electrochemical Sensors and Moisture

The OdaLog Logger Series of instruments uses an 'electro-chemical' sensor to measure levels of gas. The sensor contains an *hygroscopic* (absorbs moisture from air) electrolyte which is integral to measure gas levels. The volume of moisture inside the sensor will vary depending on the humidity levels in the air. When operating in an environment outside the sensor operating specification (outside 15-90% relative humidity) the sensor electrolyte can swell and leak in high RH, or dry out too much. High humidity is especially a concern, as a leaking sensor will result in permanent damage to the sensor. Many customers use OdaLog Loggers in enclosed 'wet-well' wastewater applications, and these environments are often above 80% humidity, and thus a special regime needs to be followed in order to ensure measurement accuracy and sensor life. The use of the OdaLog Dry Dock in this regime can greatly accelerate 'recovery time' of the instrument, reducing down-time between deployments.

SL & LL OdaLog Models

SL and LL model OdaLog Instruments use different sensors. SL sensors absorb moisture from the air quickly, and subsequently the moisture can be removed quickly. LL sensors on the other hand, absorb moisture from the air more slowly, and therefore the moisture can only be removed slowly. For highhumidity environments, the LL model is recommended, as it can be deployed for longer periods without damage (and less frequent recovery).



As a general rule of thumb, when dealing with environments which 'often' rise above 80% RH:

- * OdaLog SL models should be deployed no longer than 1 week at a time (with 3.5 days recovery on Dry Dock or 1 week in 'office' conditions)
- * OdaLog LL models should be deployed no longer than 4 weeks at a time (with 2 weeks recovery time on Dry Dock or 4 weeks in 'office' conditions)

Authorised distributor

In Australia:

For customer service, call 1300-735-292

To email an order, ordersau@thermofisher.com

To order online: thermofisher.com

In New Zealand:

For customer service, call 0800-933-966

To email an order, ordersnz@thermofisher.com

To order online: thermofisher.com

ThermoFisher
S C I E N T I F I C