

Intracellular pH (pHi) Detection

Table 1 Intracellular pH (pHi) detection products

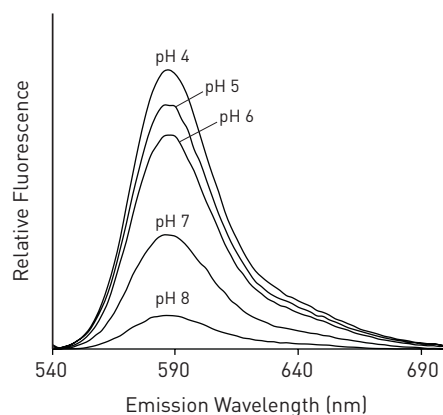
pHrodo™ Red AM Intracellular pH Indicator (Cat. no. P35372)	Amount	Concentration	Storage
pHrodo™ Red AM Intracellular pH Indicator (Component A)*	1 × 50 µL	5 mM in DMSO	<ul style="list-style-type: none"> • 2–8°C • Desiccate • Protect from light
PowerLoad™ Concentrate (Component B)	1 × 1 mL	100X	
pHrodo™ Green AM Intracellular pH Indicator (Cat. no. P35373)	Amount	Concentration	Storage
pHrodo™ Green AM Intracellular pH Indicator (Component A)*	1 × 50 µL	10 mM in DMSO	<ul style="list-style-type: none"> • 2–8°C • Desiccate • Protect from light
PowerLoad™ Concentrate (Component B)	1 × 1 mL	100X	
pHrodo™ AM Variety Pack (Cat. no. P35380)	Amount	Concentration	Storage
pHrodo™ Red AM Intracellular pH Indicator (Component A)*	1 × 10 µL	5 mM in DMSO	<ul style="list-style-type: none"> • 2–8°C • Desiccate • Protect from light
pHrodo™ Green AM Intracellular pH Indicator (Component B)*	1 × 10 µL	10 mM in DMSO	
PowerLoad™ Concentrate (Component C)	1 × 200 µL	100X	
Intracellular pH Calibration Buffer Kit (Cat. no. P35379)	Amount	Concentration	Storage
Buffer A, pH 4.5 (Component A)	1 × 50 mL	—	<ul style="list-style-type: none"> • 2–8°C • Dessiccate
Buffer B, pH 5.5 (Component B)	1 × 50 mL	—	
Buffer C, pH 6.5 (Component C)	1 × 50 mL	—	
Buffer D, pH 7.5 (Component D)	1 × 50 mL	—	
Valinomycin (Component E)	1 × 5 mg	—	
Nigericin, free acid (Component F)	1 × 5 mg	—	
DMSO, anhydrous (Component G)	1 × 1.3 mL	—	
*Approximate Fluorescence Excitation and Emission, in nm: pHrodo™ Red AM: 560/580, in nm; pHrodo™ Green AM: 509/533, in nm,			

Introduction

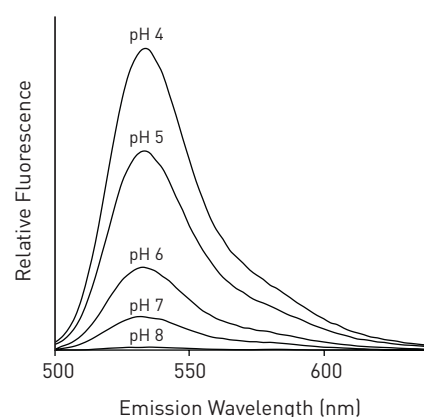
Intracellular pH (pHi) or cytosolic pH is tightly regulated within cells, and many cellular functions are dependent upon proper regulation of intracellular pH.^{1,2} Imbalance in intracellular pH can effect many cellular functions, such as ionic homeostasis, reactive oxygen species balance, apoptosis, cell cycle, and cellular mobility.³⁻⁵ pHrodo™ Red AM and pHrodo™ Green AM are novel fluorogenic probes for measuring intracellular pH in live cells. These reagents are weakly fluorescent at neutral pH, but increase their fluorescence with a drop in pH. They have pK_a of ~6.5, and are sensitive to pH change in the range of pH 9–4. Modification of pHrodo™ Red and Green dyes with AM ester groups results in an uncharged molecule that can permeate cell membranes. Once inside the cell, the lipophilic blocking groups are cleaved by nonspecific esterases, resulting in a compound that is retained within the intracellular space. With the help of Intracellular pH buffer kit, this pH change can be quantified.

- The pK_a of ~6.5 is ideally matched to the normal range of cytoplasmic pH (~6.8–7.4).
- Responses to pH across most of cellular function, in the range of pH 9–4.
- With the drop in pH, the fluorescence increases, and the reverse is true if the pH increases. This reversible property makes these dyes ideal for analyzing fluctuating cellular pH levels.
- The fluorescence emission profile of the reagents is pH-dependent (Figure 1, below), allowing the quantification of pH. Intracellular pH Kit can help in this quantification. (see **Intracellular pH Calibration**).
- Excitation and emission properties (Table 1, page 1) make these reagents useful for many platforms, including traditional fluorescence microscopy, high content screening (HCS), flow cytometry, and microplate-based fluorometry or high throughput screening (HTS).
- The acetoxymethyl (AM) ester derivative is membrane permeant, allowing noninvasive bulk loading of cells.
- Once the AM group is cleaved within the cells, the resulting compounds have excellent retention in the cells.
- These compounds are very photo stable, and their narrow spectral properties are ideal for multiplexing with other fluorophores.
- Excellent photostability and retention within the cells eliminate the need to do ratiometric measurements, allowing fluorescence reading at single excitation and emission maxima.

Figure 1 The fluorescence emission spectra of pHrodo™ Red and pHrodo™ Green conjugates

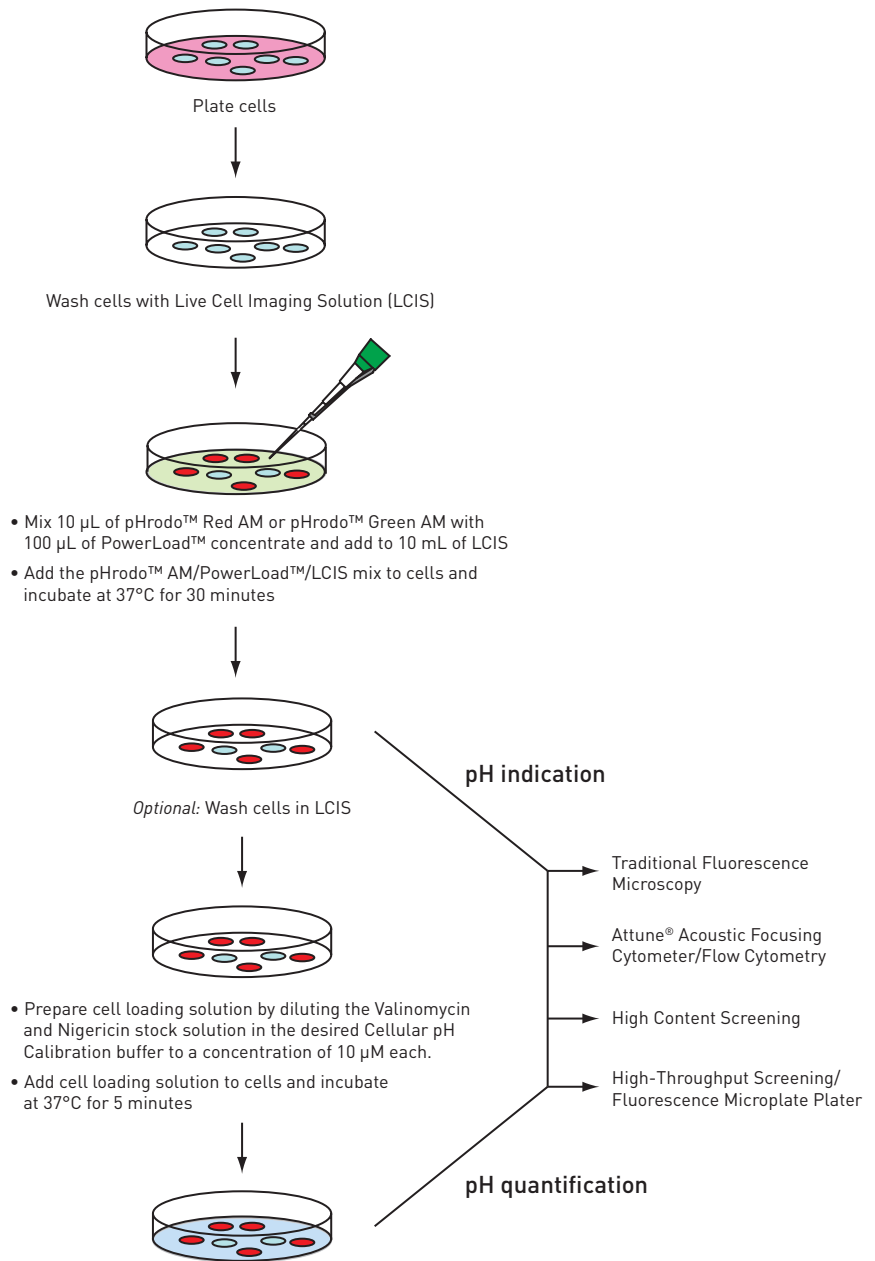


The fluorescence emission spectra of the pHrodo™ Red conjugate



The fluorescence emission spectra of the pHrodo™ Green conjugate

Figure 2 Workflow for intracellular pH (pHi) detection



pH indication: pHrodo™ Red AM and pHrodo™ Green AM

- 1.1 pHrodo™ Red AM and pHrodo™ Green AM are provided in DMSO as a 1000X solution. Thaw and mix before use.
- 1.2 Add 10 µL of the pHrodo™ dye to 100 µL of PowerLoad™ concentrate. PowerLoad™ reagent facilitates even cellular loading of AM esters; we recommend its use, but it can be omitted, if needed.
- 1.3 Dilute the dye solution into 10 mL of Live Cell Imaging Solution (LCIS) (Cat. no. A14291DJ) or other HEPES based pH 7.4 buffer to make a staining solution.
- 1.4 Remove growth medium from cells and wash once with LCIS.
- 1.5 Replace LCIS with the pHrodo™ AM staining solution.
- 1.6 Incubate at 37°C for 30 minutes.
- 1.7 Wash with LCIS if desired.
- 1.8 Analyze the cells using the appropriate instrument settings according to Ex/Em maxima in Table 1, page 1.

pH Quantification: Intracellular pH Calibration Buffer Kit

Valinomycin/Nigericin stock solution

- 2.1 Dissolve Nigericin (Component E) in 345 µL of anhydrous DMSO (Component E) to prepare a 20 mM solution.
- 2.2 Dissolve Valinomycin (Component F) in 225 µL of anhydrous DMSO (Component E) to prepare a 20 mM solution.
- 2.3 Combine 100 µL of Nigericin and 100 µL of Valinomycin solutions to prepare a 1000X Valinomycin/Nigericin stock solution (10 mM each). Aliquot and store frozen for up to 1 year.

Cellular labeling

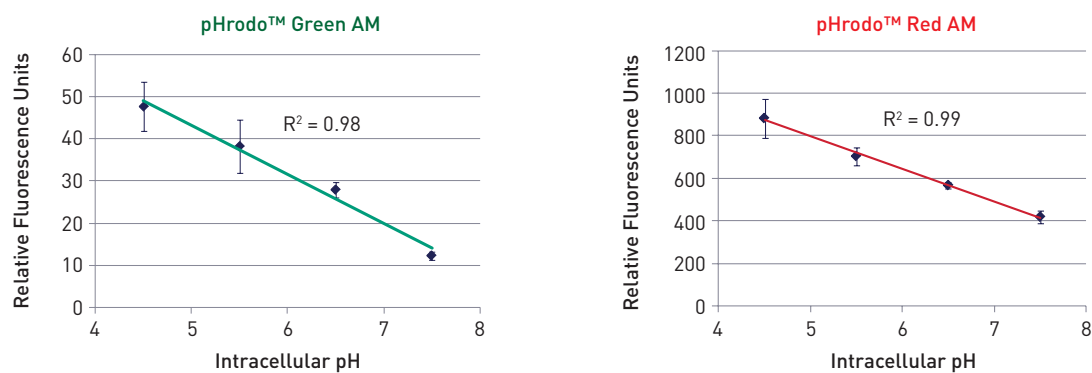
- 3.1 Prepare Cell Loading Solution by diluting 10 µL of Valinomycin/Nigericin stock solution in 10 mL of the desired Cellular pH Calibration Buffer (Component A, B, C, or D). This cell loading solution contains 10 µM of Valinomycin and 10 µM of Nigericin, in the desired pH calibration buffer.
- 3.2 Perform the desired cellular experiment with pH-sensitive pHrodo™ AM (Steps 1.1–1.5, above).
- 3.3 Wash cells 2X with Live Cell Imaging Solution (LCIS) (Cat. no. A14291DJ).
- 3.4 Replace LCIS with the 10 µM Cell Loading Solution with Valinomycin/Nigericin (from Step 3.1).
- 3.5 Incubate at 37°C for at least 5 minutes.

3.6 Analyze the cells using the appropriate Ex/Em maxima.

3.7 Repeat Steps 3.4–3.6 with the three remaining Cellular pH Calibration Buffers (Component A,B,C, or D), if desired. We recommend using at least three pH data points to get a standard curve (Figure 3, below). These data points should yield a plot, and the fluorescence value acquired in Step 3.2 can be plotted in this linear plot to get the cellular pH.

Note: Intracellular pH Calibration Buffer Kit can also be used with other conjugates of pHrodo™ and other intracellular pH indicator dyes such as SNARF® and BCECF.

Figure 3 Standard Curve using pHrodo™ Green AM and pHrodo™ Red AM with Intracellular pH Calibration Buffer Kit. U2OS cells incubated with either 10 μM pHrodo™ Green AM (left) or 5 μM pHrodo™ Red AM (right) for 30 minutes at room temperature, with PowerLoad™ reagent. Intracellular pH Calibration Buffer Kit (Cat. no. P35379) was used to clamp the intracellular pH with extracellular buffer. Four buffers included in the kit, with different pH 4.5, 5.5, 6.5, and 7.5, were used while using a high content screening (HCS) instrument. An average of three data points were plotted in the graph and a linear trendline was fitted to get the pH standard curve. Error bars represent the standard deviation.



References

1. Methods Mol Biol 637, 311 (2010);
2. Nanotechnology 24, 365 (2013);
3. Circulation 124, 1806 (2011);
4. Yonsei Med J 6, 473 (1995);
5. J Bacteriol 185, 1190 (2003).

Product List Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.	Product Name	Unit Size
P35372	pHrodo™ Red AM Intracellular pH indicator	50 µL
P35373	pHrodo™ Green AM Intracellular pH indicator	50 µL
P35379	Intracellular pH Calibration Buffer Kit	1 kit
P35380	pHrodo™ AM Variety Pack	1 kit
Related Products		
A14291DJ	Live Cell Imaging Solution	500 mL
B1150	2',7'-bis-(2-carboxyethyl)-5-(and-6)-carboxyfluorescein, acetoxymethyl ester (BCECF, AM)	1 mg
B1151	2',7'-bis-(2-carboxyethyl)-5-(and-6)-carboxyfluorescein (BCECF acid) *mixed isomers*	1 mg
B1170	2',7'-bis-(2-carboxyethyl)-5-(and-6)-carboxyfluorescein, acetoxymethyl ester (BCECF, AM) *special packaging*	20 × 50 µg
B3051	2',7'-bis-(2-carboxyethyl)-5-(and-6)-carboxyfluorescein, acetoxymethyl ester (BCECF, AM) *1 mg/mL solution in anhydrous DMSO*	1 mL
C1270	5-(and-6)-carboxy SNARF®-1	1 mg
C1271	5-(and-6)-carboxy SNARF®-1, acetoxymethyl ester, acetate	1 mg
C1272	5-(and-6)-carboxy SNARF®-1, acetoxymethyl ester, acetate *special packaging*	20 × 50 µg
C6826	5-(and-6)-chloromethyl SNARF®-1, acetate *mixed isomers* *special packaging*	20 × 50 µg
D1878	dextran, BCECF, 10,000 MW, anionic	10 mg
D1880	dextran, BCECF, 70,000 MW, anionic	10 mg
D3303	dextran, SNARF®-1, 10,000 MW, anionic	5 mg
D3304	dextran, SNARF®-1, 70,000 MW, anionic	5 mg
N1495	nigericin, free acid	10 mg
P10361	pHrodo™ Red dextran, 10,000 MW *for endocytosis*	0.5 mg
P35368	pHrodo™ Green dextran, 10,000 MW *for endocytosis*	0.5 mg

Purchaser Notification

Corporate Headquarters

5791 Van Allen Way
Carlsbad, CA 92008
USA
Phone: +1 760 603 7200
Fax: +1 760 602 6500
Email: techsupport@lifetech.com

European Headquarters

Inchinnan Business Park
3 Fountain Drive
Paisley PA4 9RF
UK
Phone: +44 141 814 6100
Toll-Free Phone: 0800 269 210
Toll-Free Tech: 0800 838 380
Fax: +44 141 814 6260
Tech Fax: +44 141 814 6117
Email: euroinfo@invitrogen.com
Email Tech: eurotech@invitrogen.com

Japanese Headquarters

LOOP-X Bldg. 6F
3-9-15, Kaigan
Minato-ku, Tokyo 108-0022
Japan
Phone: +81 3 5730 6509
Fax: +81 3 5730 6519
Email: jpinfo@invitrogen.com

Additional international offices are listed
at www.lifetechnologies.com

These high-quality reagents and materials must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Read the Safety Data Sheet provided for each product; other regulatory considerations may apply.

Obtaining Support

For the latest services and support information for all locations, go to www.lifetechnologies.com.

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support (techsupport@lifetech.com)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

SDS

Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/sds.

Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.lifetechnologies.com/support and search for the Certificate of Analysis by product lot number, which is printed on the product packaging (tube, pouch, or box).

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Disclaimer

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Important Licensing Information

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

©2013 Life Technologies Corporation. All rights reserved.