

**Dako Modified Eosin Y**
**Code CS711**
**ENGLISH**

<b>Intended use</b>	For in vitro diagnostic use. Dako Modified Eosin Y is intended for primary staining of cytoplasm (red) in formalin-fixed, paraffin-embedded tissue sections, frozen sections, and cell preparations. It is a ready-to-use reagent for use on Dako CoverStainer instrument.																																																																																																																																					
<b>Reagent provided</b>	Aqueous solution of eosin. Each bottle contains 1000 mL ready-to-use reagent. When Dako Modified Eosin Y is loaded onto Dako CoverStainer instrument, it should be used within 5 days or replaced after 3000 slides (depending on what comes first).																																																																																																																																					
<b>Precautions</b>	<ol style="list-style-type: none"> <li>1. For in vitro diagnostic use</li> <li>2. For professional users only.</li> <li>3. Wear appropriate Personal Protective Equipment to avoid contact with eyes and skin.</li> <li>4. Used and unused solution should be disposed of according to local, State and Federal regulations.</li> </ol>																																																																																																																																					
<b>Storage</b>	Store at room temperature. Do not use after expiration date stamped on bottle. If drained back into the bottle on a nightly basis, the onboard stability of the reagent is 5 days. If reagents are stored under any conditions other than those specified, the user must verify the conditions. There are no obvious signs to indicate instability of this product. If unexpected staining is observed which cannot be explained by variations in laboratory procedures and a problem with the reagent is suspected, contact Dako Technical Support.																																																																																																																																					
<b>Specimen preparation</b>	<p><u>Paraffin sections:</u> Dako Modified Eosin Y can be used for primary staining of paraffin-embedded tissue sections fixed in formalin.</p> <p><u>Frozen sections and cell preparations:</u> Dako Modified Eosin Y can be used for primary staining of acetone-fixed, frozen sections or fixed cell preparations.</p>																																																																																																																																					
<b>Staining procedure</b>	<p>Dako Modified Eosin Y, Code CS711, is ready-to-use. For use on Dako CoverStainer instrument. Please read the Administration section of the User Guide for the Dako CoverStainer instrument and program the instrument accordingly with the progressive H&amp;E staining protocol provided below.</p> <p>The recommended protocol has been validated by Dako Denmark A/S.</p> <p><b>Dako Progressive H&amp;E protocol:</b></p> <p>Pre-heating for 10 minutes at 65 °C</p> <table border="1"> <thead> <tr> <th>Step</th> <th>Dip tank</th> <th>Station</th> <th>Bottle</th> <th>Reagent</th> <th>Time min</th> <th>Time max</th> </tr> </thead> <tbody> <tr><td>1</td><td>1</td><td>1-2</td><td>A</td><td>Xylene 1*</td><td>03:00</td><td>03:30</td></tr> <tr><td>2</td><td>2</td><td>3-4</td><td>B</td><td>Xylene 2*</td><td>03:00</td><td>03:30</td></tr> <tr><td>3</td><td>3</td><td>5-6</td><td>C</td><td>96 % EtOH 1</td><td>01:00</td><td>-</td></tr> <tr><td>4</td><td>4</td><td>7-8</td><td>D</td><td>96 % EtOH 2</td><td>01:00</td><td>-</td></tr> <tr><td>5</td><td>5</td><td>9-10</td><td>E</td><td>70 % EtOH 1</td><td>02:00</td><td>-</td></tr> <tr><td>6</td><td>18</td><td>33-36</td><td>-</td><td>Tap water</td><td>01:00</td><td>-</td></tr> <tr><td>7</td><td>8</td><td>16</td><td>I</td><td>Deionized water 2</td><td>01:00</td><td>01:00</td></tr> <tr><td>8</td><td>8</td><td>15</td><td>H</td><td>Dako Harris Hematoxylin</td><td>03:00</td><td>03:00</td></tr> <tr><td>9</td><td>7</td><td>13-14</td><td>G</td><td>Deionized water 1</td><td>01:00</td><td>01:00</td></tr> <tr><td>10</td><td>9</td><td>17</td><td>J</td><td>Dako Bluing Buffer</td><td>01:00</td><td>01:00</td></tr> <tr><td>11</td><td>17</td><td>33-36</td><td>-</td><td>Tap water</td><td>01:00</td><td>01:00</td></tr> <tr><td>12</td><td>6</td><td>11-12</td><td>F</td><td>70 % EtOH 2</td><td>01:00</td><td>01:00</td></tr> <tr><td>13</td><td>10</td><td>19</td><td>L</td><td>Dako Modified Eosin Y</td><td>01:00</td><td>01:00</td></tr> <tr><td>14</td><td>11</td><td>21</td><td>N</td><td>96 % EtOH 3</td><td>01:00</td><td>01:00</td></tr> <tr><td>15</td><td>11</td><td>22</td><td>O</td><td>99.9 % EtOH 1</td><td>01:00</td><td>01:00</td></tr> <tr><td>16</td><td>12</td><td>23-24</td><td>P</td><td>99.9 % EtOH 2</td><td>01:00</td><td>01:00</td></tr> <tr><td>17</td><td>13</td><td>25-26</td><td>Q</td><td>99.9 % EtOH 3</td><td>01:00</td><td>01:00</td></tr> <tr><td>18</td><td>14-16</td><td>27-32</td><td>R, S, T</td><td>Exit solvent*</td><td>01:00</td><td>-</td></tr> </tbody> </table> <p>*Xylene or xylene substitute, Dako recommends Histo-Clear II (National Diagnostic, Inc), HS-202.</p> <p>After the staining procedure has been completed, the sections are mounted onboard Dako CoverStainer instrument using Dako CoverStainer Mounting Medium, Code CS703, or Dako Toluene-Free Mounting Medium, Code CS705. The slides must have a drying period of 10 minutes on the instrument prior to interpretation.</p> <p>It is recommended that protocol settings are verified before first use.</p>	Step	Dip tank	Station	Bottle	Reagent	Time min	Time max	1	1	1-2	A	Xylene 1*	03:00	03:30	2	2	3-4	B	Xylene 2*	03:00	03:30	3	3	5-6	C	96 % EtOH 1	01:00	-	4	4	7-8	D	96 % EtOH 2	01:00	-	5	5	9-10	E	70 % EtOH 1	02:00	-	6	18	33-36	-	Tap water	01:00	-	7	8	16	I	Deionized water 2	01:00	01:00	8	8	15	H	Dako Harris Hematoxylin	03:00	03:00	9	7	13-14	G	Deionized water 1	01:00	01:00	10	9	17	J	Dako Bluing Buffer	01:00	01:00	11	17	33-36	-	Tap water	01:00	01:00	12	6	11-12	F	70 % EtOH 2	01:00	01:00	13	10	19	L	Dako Modified Eosin Y	01:00	01:00	14	11	21	N	96 % EtOH 3	01:00	01:00	15	11	22	O	99.9 % EtOH 1	01:00	01:00	16	12	23-24	P	99.9 % EtOH 2	01:00	01:00	17	13	25-26	Q	99.9 % EtOH 3	01:00	01:00	18	14-16	27-32	R, S, T	Exit solvent*	01:00	-
Step	Dip tank	Station	Bottle	Reagent	Time min	Time max																																																																																																																																
1	1	1-2	A	Xylene 1*	03:00	03:30																																																																																																																																
2	2	3-4	B	Xylene 2*	03:00	03:30																																																																																																																																
3	3	5-6	C	96 % EtOH 1	01:00	-																																																																																																																																
4	4	7-8	D	96 % EtOH 2	01:00	-																																																																																																																																
5	5	9-10	E	70 % EtOH 1	02:00	-																																																																																																																																
6	18	33-36	-	Tap water	01:00	-																																																																																																																																
7	8	16	I	Deionized water 2	01:00	01:00																																																																																																																																
8	8	15	H	Dako Harris Hematoxylin	03:00	03:00																																																																																																																																
9	7	13-14	G	Deionized water 1	01:00	01:00																																																																																																																																
10	9	17	J	Dako Bluing Buffer	01:00	01:00																																																																																																																																
11	17	33-36	-	Tap water	01:00	01:00																																																																																																																																
12	6	11-12	F	70 % EtOH 2	01:00	01:00																																																																																																																																
13	10	19	L	Dako Modified Eosin Y	01:00	01:00																																																																																																																																
14	11	21	N	96 % EtOH 3	01:00	01:00																																																																																																																																
15	11	22	O	99.9 % EtOH 1	01:00	01:00																																																																																																																																
16	12	23-24	P	99.9 % EtOH 2	01:00	01:00																																																																																																																																
17	13	25-26	Q	99.9 % EtOH 3	01:00	01:00																																																																																																																																
18	14-16	27-32	R, S, T	Exit solvent*	01:00	-																																																																																																																																
<b>Staining interpretation</b>	Dako Modified Eosin Y stains cytoplasm of certain cells (e.g. muscle), collagen and red blood cells in various shades of pink to orange.																																																																																																																																					
<b>Additional Information</b>	Refer to User Guide for the Dako CoverStainer instrument for further information on the instrument and instructions for use.																																																																																																																																					

<b>Utilisation prévue</b>	<p>Pour utilisation diagnostique <i>in vitro</i>.</p> <p>Le réactif Dako Modified Eosin Y est destiné à la coloration primaire (en rouge) du cytoplasme sur des coupes tissulaires fixées au formol et incluses en paraffine, sur des coupes congelées et sur des préparations cellulaires. Ce réactif prêt à l'emploi est à utiliser sur l'appareil Dako CoverStainer.</p>																																																																																																																																					
<b>Réactifs fournis</b>	<p>Solution aqueuse d'éosine. Chaque flacon contient 1 000 mL de réactif prêt à l'emploi.</p> <p>Une fois chargé dans l'appareil Dako CoverStainer, le réactif Dako Modified Eosin Y doit être utilisé dans les 5 jours ou renouvelé après 3 000 lames (au premier des deux).</p>																																																																																																																																					
<b>Précautions d'emploi</b>	<ol style="list-style-type: none"> <li>1. Pour utilisation diagnostique <i>in vitro</i>.</li> <li>2. Pour utilisateurs professionnels uniquement.</li> <li>3. Porter un équipement de protection individuelle approprié pour éviter tout contact avec les yeux et la peau.</li> <li>4. Les solutions utilisées et non utilisées doivent être éliminées conformément aux réglementations locales, nationales et européennes.</li> </ol>																																																																																																																																					
<b>Conservation</b>	<p>Conserver à température ambiante. Ne pas utiliser après la date de péremption indiquée sur le flacon. S'il est remis dans son flacon chaque soir, la stabilité du réactif sur l'appareil est de 5 jours. Si les réactifs sont conservés dans des conditions autres que celles indiquées, celles-ci doivent être validées par l'utilisateur. Il n'existe pas de signe particulier pour indiquer l'instabilité de ce produit. Si une coloration inattendue est observée, qui ne peut être expliquée par des différences dans les procédures du laboratoire et qu'un problème lié au réactif est suspecté, contacter l'assistance technique de Dako.</p>																																																																																																																																					
<b>Préparation des échantillons</b>	<p><u>Coupes en paraffine :</u> Le réactif Dako Modified Eosin Y peut être utilisé pour la coloration primaire des coupes tissulaires incluses en paraffine et fixées au formol.</p> <p><u>Coupes congelées et préparations cellulaires :</u> Le réactif Dako Modified Eosin Y peut être utilisé pour la coloration primaire de coupes congelées et fixées à l'acétone ou de préparations cellulaires fixées.</p>																																																																																																																																					
<b>Procédure de coloration</b>	<p>Le réactif Dako Modified Eosin Y (code CS711) est prêt à l'emploi. Destiné à être utilisé sur l'appareil Dako CoverStainer. Lire la section Administration du guide de l'utilisateur de l'appareil Dako CoverStainer et programmer l'appareil conformément au protocole de coloration H&amp;E progressif fourni ci-dessous.</p> <p>Le protocole recommandé a été validé par Dako Denmark A/S.</p> <p><b>Protocole H&amp;E progressif de Dako :</b> Préchauffage pendant 10 minutes à 65 °C</p> <table border="1"> <thead> <tr> <th>Étape</th> <th>Réservoir</th> <th>Station</th> <th>Flacon</th> <th>Réactif</th> <th>Temps min.</th> <th>Temps max.</th> </tr> </thead> <tbody> <tr> <td>1</td> <td>1</td> <td>1-2</td> <td>A</td> <td>Xylène 1*</td> <td>03:00</td> <td>03:30</td> </tr> <tr> <td>2</td> <td>2</td> <td>3-4</td> <td>B</td> <td>Xylène 2*</td> <td>03:00</td> <td>03:30</td> </tr> <tr> <td>3</td> <td>3</td> <td>5-6</td> <td>C</td> <td>EtOH 1 à 96%</td> <td>01:00</td> <td>-</td> </tr> <tr> <td>4</td> <td>4</td> <td>7-8</td> <td>D</td> <td>EtOH 2 à 96%</td> <td>01:00</td> <td>-</td> </tr> <tr> <td>5</td> <td>5</td> <td>9-10</td> <td>E</td> <td>EtOH 1 à 70%</td> <td>02:00</td> <td>-</td> </tr> <tr> <td>6</td> <td>18</td> <td>33-36</td> <td>-</td> <td>Eau du robinet</td> <td>01:00</td> <td>-</td> </tr> <tr> <td>7</td> <td>8</td> <td>16</td> <td>I</td> <td>Eau déionisée 2</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>8</td> <td>8</td> <td>15</td> <td>H</td> <td>Dako Harris Hematoxylin</td> <td>03:00</td> <td>03:00</td> </tr> <tr> <td>9</td> <td>7</td> <td>13-14</td> <td>G</td> <td>Eau déionisée 1</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>10</td> <td>9</td> <td>17</td> <td>J</td> <td>Dako Bluing Buffer</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>11</td> <td>17</td> <td>33-36</td> <td>-</td> <td>Eau du robinet</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>12</td> <td>6</td> <td>11-12</td> <td>V</td> <td>EtOH 2 à 70%</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>13</td> <td>10</td> <td>19</td> <td>L</td> <td>Dako Modified Eosin Y</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>14</td> <td>11</td> <td>21</td> <td>N</td> <td>EtOH 3 à 96%</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>15</td> <td>11</td> <td>22</td> <td>O</td> <td>EtOH 1 à 99,9%</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>16</td> <td>12</td> <td>23-24</td> <td>P</td> <td>EtOH 2 à 99,9%</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>17</td> <td>13</td> <td>25-26</td> <td>Q</td> <td>EtOH 3 à 99,9%</td> <td>01:00</td> <td>01:00</td> </tr> <tr> <td>18</td> <td>14-16</td> <td>27-32</td> <td>R, S, T</td> <td>Solvant de sortie*</td> <td>01:00</td> <td>-</td> </tr> </tbody> </table> <p>*Xylène ou substitut de xylène, Dako recommande Histo-Clear II (National Diagnostic, Inc), HS-202.</p> <p>Une fois la procédure de coloration terminée, les sections sont installées dans l'appareil Dako CoverStainer à l'aide des milieux de montage Dako CoverStainer Mounting Medium (code CS703) ou Dako Toluene-Free Mounting Medium (code CS705). Les lames doivent sécher pendant 10 minutes sur l'instrument avant l'interprétation.</p> <p>Il est recommandé de vérifier les paramètres de protocole avant la première utilisation.</p>	Étape	Réservoir	Station	Flacon	Réactif	Temps min.	Temps max.	1	1	1-2	A	Xylène 1*	03:00	03:30	2	2	3-4	B	Xylène 2*	03:00	03:30	3	3	5-6	C	EtOH 1 à 96%	01:00	-	4	4	7-8	D	EtOH 2 à 96%	01:00	-	5	5	9-10	E	EtOH 1 à 70%	02:00	-	6	18	33-36	-	Eau du robinet	01:00	-	7	8	16	I	Eau déionisée 2	01:00	01:00	8	8	15	H	Dako Harris Hematoxylin	03:00	03:00	9	7	13-14	G	Eau déionisée 1	01:00	01:00	10	9	17	J	Dako Bluing Buffer	01:00	01:00	11	17	33-36	-	Eau du robinet	01:00	01:00	12	6	11-12	V	EtOH 2 à 70%	01:00	01:00	13	10	19	L	Dako Modified Eosin Y	01:00	01:00	14	11	21	N	EtOH 3 à 96%	01:00	01:00	15	11	22	O	EtOH 1 à 99,9%	01:00	01:00	16	12	23-24	P	EtOH 2 à 99,9%	01:00	01:00	17	13	25-26	Q	EtOH 3 à 99,9%	01:00	01:00	18	14-16	27-32	R, S, T	Solvant de sortie*	01:00	-
Étape	Réservoir	Station	Flacon	Réactif	Temps min.	Temps max.																																																																																																																																
1	1	1-2	A	Xylène 1*	03:00	03:30																																																																																																																																
2	2	3-4	B	Xylène 2*	03:00	03:30																																																																																																																																
3	3	5-6	C	EtOH 1 à 96%	01:00	-																																																																																																																																
4	4	7-8	D	EtOH 2 à 96%	01:00	-																																																																																																																																
5	5	9-10	E	EtOH 1 à 70%	02:00	-																																																																																																																																
6	18	33-36	-	Eau du robinet	01:00	-																																																																																																																																
7	8	16	I	Eau déionisée 2	01:00	01:00																																																																																																																																
8	8	15	H	Dako Harris Hematoxylin	03:00	03:00																																																																																																																																
9	7	13-14	G	Eau déionisée 1	01:00	01:00																																																																																																																																
10	9	17	J	Dako Bluing Buffer	01:00	01:00																																																																																																																																
11	17	33-36	-	Eau du robinet	01:00	01:00																																																																																																																																
12	6	11-12	V	EtOH 2 à 70%	01:00	01:00																																																																																																																																
13	10	19	L	Dako Modified Eosin Y	01:00	01:00																																																																																																																																
14	11	21	N	EtOH 3 à 96%	01:00	01:00																																																																																																																																
15	11	22	O	EtOH 1 à 99,9%	01:00	01:00																																																																																																																																
16	12	23-24	P	EtOH 2 à 99,9%	01:00	01:00																																																																																																																																
17	13	25-26	Q	EtOH 3 à 99,9%	01:00	01:00																																																																																																																																
18	14-16	27-32	R, S, T	Solvant de sortie*	01:00	-																																																																																																																																
<b>Interprétation de la coloration</b>	<p>Le réactif Dako Modified Eosin Y colore le cytoplasme de certaines cellules (par ex., du muscle), le collagène et les hématies en des teintes variées allant du rose à l'orangé.</p>																																																																																																																																					
<b>Informations supplémentaires</b>	<p>Consulter le guide de l'utilisateur de l'appareil Dako CoverStainer pour des informations supplémentaires sur l'appareil et le mode d'emploi.</p>																																																																																																																																					

- Verwendungszweck** Zur In-vitro-Diagnostik.  
Dako Modified Eosin Y ist zur primären Färbung von Zytoplasma (rot) in formalinfixierten, paraffineingebetteten Gewebeschnitten, Gefriergewebeschnitten und Zellpräparaten bestimmt. Es handelt sich um ein gebrauchsfertiges Reagenz für die Verwendung mit dem Dako CoverStainer.
- Geliefertes Reagenz** Wässrige Eosinlösung. Jede Flasche enthält 1000 mL gebrauchsfertiges Reagenz.  
Dako Modified Eosin Y ist nach dem Beladen in den Dako CoverStainer nach 5 Tagen bzw. nach 3.000 Objektträgern auszutauschen (ja nachdem, was zuerst eintritt).
- Vorsichtsmaßnahmen**
1. Zur In-vitro-Diagnostik.
  2. Nur für Fachpersonal bestimmt.
  3. Entsprechende persönliche Schutzausrüstung (PSA) tragen, um Augen- und Hautkontakt zu vermeiden.
  4. Verwendete und auch nicht verwendete Lösung ist entsprechend örtlichen, staatlichen und EU-rechtlichen Richtlinien zu entsorgen.
- Lagerung** Bei Raumtemperatur aufbewahren. Nach Ablauf des auf der Flasche aufgedruckten Verfallsdatums nicht mehr verwenden. Beim Zurückfüllen des Reagenz in die Flasche zur Aufbewahrung über Nacht beträgt die Stabilität im Gerät 5 Tage. Sollten die Reagenzien unter anderen Bedingungen als den beschriebenen aufbewahrt worden sein, so müssen diese vom Anwender verifiziert werden. Es gibt keine offensichtlichen Anhaltspunkte für die mögliche Instabilität dieses Produkts. Wenn eine unerwartete Anfärbung beobachtet wird, die durch Abweichungen in den Labormethoden nicht erklärt werden kann, und falls Verdacht auf ein Problem mit dem Reagenz besteht, bitte Kontakt mit dem technischen Kundendienst von Dako aufnehmen.
- Gewebevorbereitung**
- Paraffinschnitte:  
Dako Modified Eosin Y kann für die primäre Färbung von formalinfixierten, paraffineingebetteten Gewebeschnitten verwendet werden.
- Gefrierschnitte und Zellpräparate:  
Dako Modified Eosin Y Hematoxylin kann zur primären Färbung von azetonfixierten Gefrierschnitten oder fixierten Zellausstrichen verwendet werden.
- Färbeverfahren** Dako Modified Eosin Y, Code-Nr. CS711, ist gebrauchsfertig. Für die Verwendung mit dem Dako CoverStainer. Bitte im Benutzerhandbuch Dako CoverStainer, Abschnitt „Administration“ nachschlagen, und das Gerät auf das unten aufgeführte progressive H&E-Färbeprotokoll programmieren.  
Das empfohlene Protokoll wurde von Dako Denmark A/S getestet.

**Progressives H&E-Protokoll von Dako**

Vorwärmen für 10 Minuten bei 65 °C

Schritt	Eintauchtank	Station	Flasche	Reagenz	Min.-Zeit	Max.-Zeit
1	1	1-2	A	Xylol 1*	03:00	03:30
2	2	3-4	B	Xylol 2*	03:00	03:30
3	3	5-6	C	96% EtOH 1	01:00	–
4	4	7-8	D	96% EtOH 2	01:00	–
5	5	9-10	E	70% EtOH 1	02:00	–
6	18	33-36	–	Leitungswasser	01:00	–
7	8	16	I	Entionisiertes Wasser 2	01:00	01:00
8	8	15	H	Dako Harris Hematoxylin	03:00	03:00
9	7	13-14	G	Entionisiertes Wasser 1	01:00	01:00
10	9	17	J	Dako Bläuungspuffer	01:00	01:00
11	17	33-36	–	Leitungswasser	01:00	01:00
12	6	11-12	F	70% EtOH 2	01:00	01:00
13	10	19	L	Dako Modified Eosin Y	01:00	01:00
14	11	21	n	96% EtOH 3	01:00	01:00
15	11	22	O	99,9% EtOH 1	01:00	01:00
16	12	23-24	P	99,9% EtOH 2	01:00	01:00
17	13	25-26	Q	99,9% EtOH 3	01:00	01:00
18	14-16	27-32	R, S, T	Ausgabelösungsmittel*	01:00	–

\*Xylol oder Xylolersatzstoff, Dako empfiehlt Histo-Clear II (National Diagnostic, Inc), HS-202.

Nach Abschluss des Färbeverfahrens werden die Schnitte mithilfe von Dako CoverStainer Mounting Medium, Code-Nr. CS703, oder Dako Toluene-Free Mounting Medium, Code-Nr. CS705, in den Dako CoverStainer beladen. Vor einer Auswertung müssen die Objektträger 10 Minuten lang auf dem Instrument trocknen.

Es wird empfohlen, vor der erstmaligen Verwendung die Protokolleinstellungen zu überprüfen.

- Auswertung der Färbung** Dako Modified Eosin Y färbt Zytoplasma bestimmter Zellen (z. B. Muskelzellen), Kollagen und Erythrozyten in verschiedenen Farbtönen von Pink bis Orange.
- Weitere Informationen** Weitere Informationen über das Gerät und die Gebrauchsanweisung finden Sie im Benutzerhandbuch für das Dako CoverStainer-Gerät.

---

**Companion Products/ Produits d'accompagnement/ Begleitprodukte**


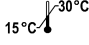





Code CS709, Dako Harris Hematoxylin  
Code CS702, Dako Bluing Buffer

Code CS703, Dako Mounting Medium  
or/ ou/ oder  
Code CS705, Dako Toluene Free Mounting Medium

Code CS704, Dako Cover Glass

---

**Explanation of symbols/ Explications des symboles/ Erläuterung der Symbole**

 Catalogue number Référence du catalogue Katalognummer	 15°C - 30°C Temperature limitation Limites de température Zulässiger Temperaturbereich	 Manufacturer Fabricant Hersteller
 In vitro diagnostic medical device Dispositif médical de diagnostic in vitro In-vitro-Diagnostikum	 Batch code Code du lot Chargenbezeichnung	
 Consult instructions for use Consulter le mode d'emploi Gebrauchsanweisung beachten	 Use by Utiliser avant Verwendbar bis	

Manufactured by: / Fabriqué par : / Hergestellt von:  
American Master Tech Scientific (AMTS), Inc, USA

Manufactured for: / Fabriqué pour : / Hergestellt für:  
Dako Denmark A/S, An Agilent Technologies Company