

References: Influenza Neuraminidase Activity Quantitation with NA-Star® Chemiluminescent Substrate/NA-Star® Influenza Neuraminidase Inhibitor Resistance Detection Kit

- Buxton, RC, B Edwards, RR Juo, JC Voyta, M Tisdale and RC Bethell (2000). Development of a sensitive chemiluminescent neuraminidase assay for the determination of influenza virus susceptibility to Zanamivir. *Anal Biochem* 280:291-300. doi:10.1006/abio.2000.4517.
- Boivin, G and N Goyette (2002). Susceptibility of recent Canadian influenza A and B virus isolates to different neuraminidase inhibitors. *Antiviral Research* 54:143-147.
- Aymard, M, O Ferraris, L Gerentes, J Jolly and N Kessler (2003). Neuraminidase assays. *Dev Biol* 115:75-83.
- McKimm-Breschkin, J, T Trivedi, A Hampson, A Hay, A Klimov, M Tashiro, F Hayden and M Zambon (2003). Neuraminidase sequence analysis and susceptibilities of influenza virus clinical isolates to Zanamivir and Oseltamivir. *Antimicrobial Agents and Chemotherapy* 47(7):2264-2272. doi:10.1128/AAC.47.7.2264-2272.2003.
- Mungall, BA, X Xu and A Klimov (2003). Assaying susceptibility of avian and other influenza A viruses to Zanamivir: Comparisons of fluorescent and chemiluminescent neuraminidase assays. *Avian Diseases* 47:1141-1144.
- Wetherall, NT, T Trivedi, J Zeller, C Hodges-Savola, JL McKimm-Breschkin, M Zambon and FG Hayden (2003). Evaluation of neuraminidase enzyme assays using different substrates to measure susceptibility of influenza virus clinical isolates to neuraminidase inhibitors: Report of the Neuraminidase Inhibitor Susceptibility Network. *J Clin Microbiol* 41(2):742-750. doi:10.1128/JCM.41.2.742-750.2003.
- Hurt, AC, JL McKimm-Breschkin, M McDonald, IG Barr, N Komadina and AW Hampson (2004). Identification of a human influenza type B strain with reduced sensitivity to neuraminidase inhibitor drugs. *Virus Research* 103:205-211. doi:10.1016/j.virusres.2004.02.035.
- Mungall BA, X Xu and A Klimov (2004). Surveillance of influenza isolates for susceptibility to neuraminidase inhibitors during the 2000-2002 influenza seasons. *Virus Research* 103:195-197. doi:10.1016/j.virusres.2004.02.033.
- Jackson, D, W Barclay and T Zurcher (2005). Characterization of recombinant influenza B viruses with key neuraminidase inhibitor resistance mutations. *J Antimicrobial Chemotherapy* 55(2):162-169. doi:10.1093/jac/dkh528.
- The World Health Organization Global Influenza Program Surveillance Network (2005). Evolution of H5N1 avian influenza viruses in Asia. *Emerging Infectious Diseases* 11(10):1515-1521.
- Monto, AS, JL McKimm-Breschkin, C Macken, AW Hampson, A Hay, A Klimov, M Tashiro, RG Webster, M Aymard, FG Hayden and M Zambon (2006). Detection of influenza viruses resistant to neuraminidase inhibitors in global surveillance during the first 3 years of their use. *Antimicrobial Agents and Chemotherapy* 50(7):2395-2402. doi:10.1128/AAC.01339-05.
- Zurcher, T, PJ Yates, J Daly, A Sahasrabudhe, M Walters, L Dash, M Tisdale and JL McKimm-Breschkin (2006). Mutations conferring zanamivir resistance in human influenza virus N2 neuraminidases compromise virus fitness and are not stably maintained *in vitro*. *J Antimicrobial Chemotherapy* 58:723-732. doi:10.1093/jac/dk1321.
- Bauer, K, C Schrader, J Suess, P Wutzler and M Schmidtke (2007). Neuraminidase inhibitor susceptibility of porcine H3N2 influenza A viruses isolated in Germany between 1982 and 1999. *Antiviral Res* 75:219-226. doi:10.1016/j.antiviral.2007.03.007.
- Bright, RA, DM Carter, S Daniluk, FR Toapanta, A Ahmad, V Gavrilov, M Massare, P Pushko, N Mytle, T Rowe, G Smith and TM Ross (2007). Influenza virus-like particles elicit broader immune responses than whole virion inactivated influenza virus or recombinant hemagglutinin. *Vaccine* 25:3871-3878. doi:10.1016/j.vaccine.2007.01.106.
- Bouvier, NM, AC Lowen and P Palese (2008). Oseltamivir-resistant influenza A viruses are transmitted efficiently among guinea pigs by direct contact but not by aerosol. *J Virol* 82(20):10052-10058. doi:10.1128/JVI.01226-08.
- Le, MTQ, HFL Wertheim, HD Nguyen, W Taylor, PVM Hoang, et al. (2008). Influenza A H5N1 clade 2.3.4 virus with a different antiviral susceptibility profile replaced clade 1 virus in humans in northern Vietnam. *PloS ONE* 3(10):e3339. doi:10.1371/journal.pone.0003339.
- Sheu, TG, VM Deyde, M Okomo-Adhiambo, RJ Garten, X Xu, RA Bright, EN Butler, TR Wallis, AI Klimov and LV Gubareva (2008). Surveillance for neuraminidase inhibitor resistance among human influenza A and B viruses circulating worldwide from 2004 to 2008. *Antimicrobial Agents Chemotherapy* 52(9):3284-3292. doi:10.1128/AAC.00555-08.

- Taylor, WRJ, BN Thinh, GT Anh, P Horby, H Wertheim, et al. (2008). Oseltamivir is adequately absorbed following nasogastric administration to adult patients with severe H5N1 influenza. *PLoS ONE* 3(10): e3410. doi:10.1371/journal.pone.0003410.
- Aljofan, M, ML Sganga, MK Lo, CL Rootes, M Porotto, AG Meyer, S Saubern, A Moscona and BA Mungall (2009). Antiviral activity of gliotoxin, gentian violet and brilliant green against Nipah and Hendra virus *in vitro*. *Virology J* 6:187. doi:10.1186/1743-422X-6-187.
- An, J, DCW Lee, AHY Law, CLH Yang, LLM Poon, ASY Lau and SJM Jones (2009). A novel small-molecule inhibitor of the avian influenza H5N1 virus determined through computational screening against the neuraminidase. *J Med Chem* 52:2667-2672.
- Bauer, K, M Richter, P Wutzler and M Schmidtke (2009). Different neuraminidase inhibitor susceptibilities of human H1N1, H1N2, and H3N2 influenza A viruses isolated in Germany from 2001 to 2005/2006. *Antiviral Res* 82:34-41. doi:10.1016/j.antiviral.2009.01.006.
- Chen, H, CL Cheung, H Tai, P Zhao, JFW Chan, VCC Cheng, K-H Chan and K-Y Yuen (2009). Oseltamivir-resistant influenza A pandemic (H1N1) 2009 virus, Hong Kong, China. *Emerg Infect Dis* 15(12):1970-1972. doi:10.3201/eid1512.091057.
- Cheng, PKC, TWC Leung, ECM Ho, PCK Leung, AYY Ng, MYY Lai and WWL Lin (2009). Oseltamivir- and amantadine-resistant influenza viruses A (H1N1). *Emerg Infect Dis* 15(6):966-968. doi:10.3201/eid1506.081357.
- Deyde, VM, M Okomo-Adhiambo, TG Sheu, TR Wallis, A Fry, N Dharan, AI Klimov and LV Gubareva (2009). Pyrosequencing as a tool to detect molecular markers of resistance to neuraminidase inhibitors in seasonal influenza A viruses. *Antiviral Res* 81:16-24. doi:10.1016/j.antiviral.2008.08.008.
- Dharan, NJ, LV Gubareva, JJ Meyer, M Okomo-Adhiambo, RC McClinton, SA Marshall, K St George, S Epperson, L Brammer, AI Klimov, JS Bresee and AM Fry (for the Oseltamivir-Resistance Working Group) (2009). Infections with oseltamivir-resistant influenza A (H1N1) virus in the United States. *J Am Med Assoc* 301(10), doi:10.1001/jama.2009.294.
- Hashem, AM, AS Flaman, A Farnsworth, EG Brown, G Van Domselaar, R He and X Li (2009). Aurintricarboxylic acid is a potent inhibitor of influenza A and B virus neuraminidases. *PLoS ONE* 4(12): e8350. doi:10.1371/journal.pone.0008350.
- Hung, H-C, C-P Tseng, J-M Yang, Y-W Ju, S-N Tseng, Y-F Chen, Y-S Chao, H-P Hsieh, S-R Shih and JT-A Hsu (2009). Aurintricarboxylic acid inhibits influenza virus neuraminidase. *Antiviral Res* 81:123-131. doi:10.1016/j.antiviral.2008.10.006.
- Kawakami, C, M Obuchi, M Saikusa, Y Noguchi, M Ujike, T Odagiri, M Iwata, T Toyozawa and M Tashiro (2009). Isolation of oseltamivir-resistant influenza A/H1N1 virus of different origins in Yokohama City, Japan, during the 2007-2008 influenza season. *Jpn J Infect Dis* 62:83-86.
- Leung, TWC, ALS Tai, PKC Cheung, MSY Kong and W Lim (2009). Detection of an oseltamivir-resistant pandemic influenza A/H1N1 virus in Hong Kong. *J Clin Virol* 46:298-299. doi:10.1016/j.jcv.2009.08.004.
- Lin, X, J Zhou, Y Zhang, J Wu, F Zhang, Z Li, Y Zhang, S Bi, Y Shu and Y Wang (2009). Oseltamivir boosts 2009 H1N1 virus infectivity *in vitro*. *Biochem Biophys Res Comm* 390:1305-1308. doi:10.1016/j.bbrc.2009.10.142.
- Novel Swine-Origin Influenza A (H1N1) Virus Investigation Team (2009). Emergence of a novel swine-origin influenza A (H1N1) virus in humans. *N Engl J Med* 360:2605-2615. doi:10.106/NEJMoa0903810.
- Shinde, V, et al. (2009) Triple-reassortant swine influenza A (H1) in humans in the United States, 2005-2009. *N Engl J Med* 360:2616-2625. doi:10.1056/NEJMoa0903812.
- Tashiro, M, JL McKimm-Breschkin, T Saito, A Klimov, C Macken, M Zambon and FG Hayden for NISN (2009). Surveillance for neuraminidase-inhibitor-resistant influenza viruses in Japan, 1996-2007. *Antiviral Therapy* 14:751-761. doi:10.3851/IMP1194.
- Triana-Baltzer, GB, LV Gubareva, AI Klimov, DF Wurtman, RB Moss, M Hedlund, JL Larson, RB Belshe and F Fang (2009). Inhibition of neuraminidase inhibitor-resistant influenza virus by DAS181, a novel sialidase fusion protein. *PLoS ONE* 4(11): e7838. doi:10.1371/journal.pone.0007838.
- Wen, W-H, M Lin, C-Y Su, S-Y Wang, Y-SE Cheng, J-M Fang and C-H Wong (2009). Synergistic effect of Zanamivir-porphyrin conjugates on inhibition of neuraminidase and inactivation of influenza virus. *J Med Chem* 52(15):4903-4910. doi:10.1021/jm900515g.
- Bastien, N, NA Antonishyn, K Brandt, CE Wong, K Chokani, N Vegh, GB Horsman, S Tyler, MR Graham, FA Plummer, PN Levett and Y Li (2010). Human infection with a triple-reassortment swine influenza A(H1N1) virus containing the hemagglutinin and neuraminidase genes of seasonal influenza virus. *J Infectious Diseases* 201:1178-1182. doi:10.1086/651507.

- Blair, PJ, TF Wierzba, S Touch, S Vonthanak, X. Xu, RJ Garten, MA Okomo-Adhiambo, AI Klimov, MR Kasper and SD Putnam (2010). Influenza epidemiology and characterization of influenza viruses in patients seeking treatment for acute fever in Cambodia. *Epidemiol Infect* 138:199-209. doi:10.1017/S095026880999063X.
- Grienke, U, M Schmidtke, J Kirchmair, K Pfarr, P Wutzler, R Durrwald, G Wolber, KR Liedl, H Stuppner and JM Rollinger (2010). Antiviral potential and molecular insight into neuraminidase inhibiting diarylheptanoids from *Alpinia katsumadai*. *J Med Chem* 53:778-786. doi:10.1021/jm901440f.
- Hassantoufighi, A, H Zhang, M Sandbulte, J Gao, J Manischewitz, L King, H Golding, TM Straight and MC Eichelberger (2010). A practical influenza neutralization assay to simultaneously quantify hemagglutinin and neuraminidase-inhibiting antibody responses. *Vaccine* 28:790-797. doi:10.1016/j.vaccine.2009.10.066.
- Jonges, M, WM Liu, E van der Vries, R Jacobi, I Pronk, C Boog, M Koopmans, A Meijer and E Soethout (2010). Influenza virus inactivation for studies of antigenicity and phenotypic neuraminidase inhibitor resistance profiling. *J Clin Microbiol* 48(3):928-940. doi:10.1128/JCM.02045-09.
- Lai, JCC, WWL Chan, F Kien, JM Nicholls, JSM Peiris and J-M Garcia (2010). Formation of virus-like particles from human cell lines exclusively expressing influenza neuraminidase. *J General Virology* 91:2322-2330. doi:10.1099/vir.0.019935-0.
- Matsuzaki, Y, K Mizuta, Y Aoki, A Suto, C Abiko, K Sanjoh, K Sugawara, E Takashita, T Itagaki, Y Katsushima, M Ujike, M Obuschi, T Odagiri and M Tashiro (2010). A two-year survey of the oseltamivir-resistant influenza A(H1N1) virus in Yamagata, Japan and clinical effectiveness of oseltamivir and zanamivir. *Virology Journal* 7:53. <http://www.virologyj.com/content/7/1/53>.
- Nguyen, HT, TG Sheu, VP Mishin, AI Klimov and LV Gubareva (2010). Assessment of pandemic and seasonal influenza A (H1N1) virus susceptibility to neuraminidase inhibitors in three enzyme activity inhibition assays. *Antimicrobial Agents Chemotherapy* 54(9):3671-3677. doi:10.1128/AAC.00581-10.
- Okomo-Adhiambo, M, K Sleeman, K Ballenger, HT Nguyen, VP Mishin, TG Sheu, J Smagala, Y Li, AI Klimov and LV Gubareva (2010). Neuraminidase inhibitor susceptibility testing in human influenza viruses: a laboratory surveillance perspective. *Viruses* 2:2269-2289. doi:10.3390/v2102269.
- Seibert, CW, M Kaminski, J Philipp, D Rubbenstroth, RA Albrecht, F Schwalm, S Stertz, RA Medina, G Kochs, A Garcia-Sastre, P Staeheli and P Palese (2010). Oseltamivir-resistant variants of the 2009 pandemic H1N1 influenza A virus are not attenuated in the guinea pig and ferret transmission models. *J Virology* 84(21):11219-11226. doi:10.1128/JVI.01424-10.
- Smee, DF, BL Hurst, M-H Wong, KW Bailey, EB Tarbet, JD Morrey and Y Furuta (2010). Effects of the combination of favipiravir (T-705) and oseltamivir on influenza A virus infections in mice. *Antimicrobial Agents and Chemotherapy* 54(1):126-133. doi:10.1128/AAC.00933-09.
- Yang, J-R, Y-P Huang, Y-Cheng Lin, C-H Su, C-Y Kuo, L-C Hsu, H-S Wu and M-T Liu (2010). Early findings of oseltamivir-resistant pandemic (H1N1) 2009 influenza A viruses in Taiwan. *Antiviral Res*, in press. doi:10.1016/j.antiviral.2010.09.014.
- Zhang, Y, X Lin, G Wang, J Zhou, J Lu et al. (2010). Neuraminidase and hemagglutinin matching patterns of a highly pathogenic avian and two pandemic H1N1 influenza A viruses. *PLoS ONE* 5(2): e9167. doi:10.1371/journal.pone.0009167.

For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.

© 2010 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.