

Selected Literature References:

- Detection of free radicals in beer oxidation: H. Kaneda, et al.; J. Food Sci. **53**, 885 (1988)
- The role of free radicals in beer oxidation: H. Kaneda, et al.; J. Am. Soc. Brew. Chem. **47**, 49 (1989)
- Improvement for oxidative flavor stability of beer - role of OH-radical in beer oxidation: M. Uchida & M. Ono; J. Am. Soc. Brew. Chem. **54**, 198 (1996)
- Electron spin resonance spin trapping identification of radicals formed during aerobic forced aging of beer: M.L. Andersen & L.H. Skibsted; J. Agric. Food Chem. **46**, 1272 (1998)
- The role of riboflavin in beer flavour instability : EPR studies and the application of flavin binding proteins: C. Laane, et al.; J. Inst. Brew. **105**, 392 (1999)
- The stability of beer characterized by means of radical traps: A. Stasko, et al.; Kvasny Prumysl **45**, 315 (1999)
- Characterization of beer stability by means of spin trapping (an EPR study): A. Stasko, et al.; Monatsschr. Brauwiss. **53**, 4 (2000)
- Potential antioxidants in beer assessed by ESR spin trapping: M.L. Andersen, et al.; J. Agric. Food Chem. **48**, 3106 (2000)
- Modification of the levels of polyphenols in wort and beer by addition of hexamethylenetetramine or sulfite during mashing: M.L. Andersen & L.H. Skibsted; J. Agric. Food Chem. **49**, 5232 (2001)
- Electron paramagnetic resonance (EPR) profiling for potential flavor stability improvements in beer: R.T. Foster, et al.; MBAA TQ **38**, 247 (2001)
- The influence of additives on beer stability investigated by EPR spectroscopy: V. Brezova, et al.; Spectrochimica Acta (Part A) **58**, 1279 (2002)
- Fundamentals on electro spin resonance spectroscopy (ESR) and research about the correlation between oxidative beer stability and sulphite content: T. Kunz, et al.; Monatsschr. Brauwiss. **55**, 140 (2002)
- Determination of antioxidants in brewing: some aspects about the use of selected chemical and physical assays: A. Stephan, et al.; Eur. Brew. Conv. **31**, 99 (2002)
- Radicaloid-type oxidative decomposition of beer bittering agents revealed: K. Huvaere, et al.; Chemistry **9**, 4693 (2003)
- Stability index: A new approach to measure the flavor stability of beer: O. Franz & W. Back; MBAA TQ **40**, 20 (2003)
- Photooxidative degradation of beer bittering principles: a key step on the route to lightstruck flavor formation in beer: K. Huvaere, et al.; J. Agric. Food Chem. **53**, 1489 (2005)
- Antioxidant pool in beer and kinetics of EPR spin-trapping: N.M. Kocherginsky, et al.; J. Agric. Food Chem. **53**, 6870 (2005)